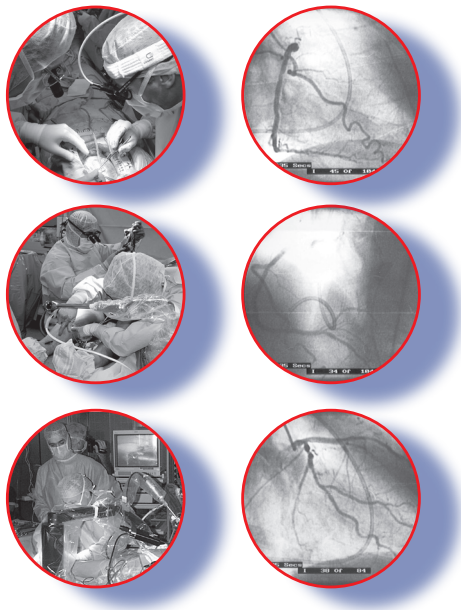




# ADMA-ELISA



**The classic test and  
gold standard for  
more than 10 years**

**List of published studies using the  
ADMA-ELISA of DLD Diagnostika GmbH**



## 2014

last update June 2014

- Beltrán, LM; Muñoz Hernández, R; de Pablo Bernal, RS; García Morillo, JS; Egido, J; Noval, ML; Ferrando-Martinez, S; Blanco-Colio, LM; Genebat, M; Villar, JR; Moreno-Luna, R; Moreno, JA  
**Reduced sTWEAK and increased sCD163 levels in HIV-infected patients: modulation by antiretroviral treatment, HIV replication and HCV co-infection**  
PLoS ONE 2014; 9: e90541
  - Borg, MK; Ivarsen, P; Brøndum, E; Povlsen, JV; Aalkjær, C  
**Hemodynamics and function of resistance arteries in healthy persons and end stage renal disease patients**  
PLoS ONE 2014; 9: e94638
  - Chiavaroli, V; Desses, L; Giorgis, TD; Giannini, C; Marcovecchio, ML; Chiarelli, F; Mohn, A  
**Is Asymmetric Dimethylarginine Associated with Being Born Small and Large for Gestational Age?**  
Antioxid. Redox Signal. 2014; :
  - Dayem, SM; Battah, AA; El-Shehaby, A; Bohy, AE  
**Asymmetric dimethyl L-arginine, nitric oxide and cardiovascular disease in adolescent type 1 diabetics**  
J. Pediatr. Endocrinol. Metab. 2014; :
  - Deftereos, S; Bouras, G; Tsounis, D; Papadimitriou, C; Hatzis, G; Raisakis, K; Panagopoulou, V; Kaoukis, A; Ioannidis, A; Deftereos, G; Kossyvakis, C; Manolis, AS; Alexopoulos, D; Stefanadis, C; Cleman, MW; Giannopoulos, G  
**Association of asymmetric dimethylarginine levels with treadmill-stress-test-derived prognosticators**  
Clin. Biochem. 2014; :
  - Ghebremariam, YT; Erlanson, DA; Cooke, JP  
**A novel and potent inhibitor of dimethylarginine dimethylaminohydrolase: a modulator of cardiovascular nitric oxide**  
J. Pharmacol. Exp. Ther. 2014; 348: 69-76
  - Karakurt, F; Carlioglu, A; Kaygusuz, I; Gumus, II; Uz, B; Akdeniz, D;  
**Effect of ethinyl estradiol-cyproterone acetate treatment on asymmetric dimethyl-arginine levels in women with polycystic ovary syndrome**  
Arch. Gynecol. Obstet. 2014; 289: 135-140
  - Petr, J; Michal, H; Jan, S; Ladislav, T; Jiri, K; Josef, S  
**Reactive hyperaemia index as a marker of endothelial dysfunction in children with Crohn's disease is significantly lower than healthy controls**  
Acta Paediatr. 2014; 103: 55-60
  - Začiragić, A; Huskić, J; Mulabegović, N; Avdagić, N; Valjevac, A; Hasić, S; Jadrić, R;  
**An assessment of correlation between serum asymmetric dimethylarginine and glycated haemoglobin in patients with type 2 diabetes mellitus**  
Bosn J Basic Med Sci 2014; 14:21-24
- ## 2013
- Bakr, A; Pak, O; Taye, A; Hamada, F; Hemeida, R; Janssen, W; Gierhardt, M; Ghofrani, HA; Seeger, W; Grimminger, F; Schermuly, RT; Witzentrath, M; Brandes, RP; Huang, N; Cooke, JP; Weissmann, N; Sommer, N  
**Effects of dimethylarginine dimethylaminohydrolase-1 overexpression on the response of the pulmonary vasculature to hypoxia**  
Am. J. Respir. Cell Mol. Biol. 2013; 49: 491-500
  - Betz, B; Möller-Ehrlich, K; Kress, T; Kniepert, J; Schwedhelm, E; Böger, RH; Wanner, C; Sauvant, C; Schneider, R  
**Increased symmetrical dimethylarginine in ischemic acute kidney injury as a causative factor of renal L-arginine deficiency**  
Transl Res 2013; 162: 67-76
  - Damasceno, NR; Sala-Vila, A; Cofán, M; Pérez-Heras, AM; Fitó, M; Ruiz-Gutiérrez, V; Martínez-González, MÁ; Corella, D; Arós, F; Estruch, R; Ros, E  
**Mediterranean diet supplemented with nuts reduces waist circumference and shifts lipoprotein subfractions to a less atherogenic pattern in subjects at high cardiovascular risk**  
Atherosclerosis 2013; 230: 347-353
  - Eisinger, K; Krautbauer, S; Wiest, R; Karrasch, T; Hader, Y; Scherer, MN; Farkas, S; Aslanidis, C; Buechler, C  
**Portal vein omentin is increased in patients with liver cirrhosis but is not associated with complications of portal hypertension**  
Eur. J. Clin. Invest. 2013; 43: 926-932
  - Eissa, SS  
**Serum Asymmetric Dimethylarginine (ADMA) and Insulin Resistance in Polycystic Ovary Syndrome**  
Journal of Applied Sciences Research 2013; 9:
  - Eleuterio, NM; Palei, AC; Rangel Machado, JS; Tanus-Santos, JE; Cavalli, RC; Sandrim, VC  
**Relationship between adiponectin and nitrite in healthy and preeclampsia pregnancies**  
Clin. Chim. Acta 2013; 423: 112-115
  - Eloot, S; Van Biesen, W; Glorieux, G; Neiryck, N; Dhondt, A; Vanholder, R  
**Does the adequacy parameter Kt/V(urea) reflect uremic toxin concentrations in hemodialysis patients?**  
PLoS ONE 2013; 8: e76838

- Esfahani, PS; Gharakhanlou, R; Karimian, J; Khazaei, M; Feizi, A; Safarzade, A  
**Effect of resistance training on plasma nitric oxide and asymmetric dimethylarginine concentrations in type I diabetic rats**  
Int J Prev Med 2013; 4: S78-84
- Ghebremariam, YT; Yamada, K; Lee, JC; Johnson, CL; Atzler, D; Anderssohn, M; Agrawal, R; Higgins, JP; Patterson, AJ; Böger, RH; Cooke, JP  
**FXR agonist INT-747 upregulates DDAH expression and enhances insulin sensitivity in high-salt fed Dahl rats**  
PLoS ONE 2013; 8: e60653
- Gürel, E; Tigen, K; Karaahmet, T; Geçmen, C; Mutlu, B; Başaran, Y  
**Predictive value of plasma asymmetric dimethylarginine, homocysteine, and high-sensitive CRP levels in occult coronary artery disease : A multidetector-row computed tomography study**  
Herz 2013; null
- Hudson, CL; Zemlin, AE; Ipp, H  
**The cardiovascular risk marker asymmetric dimethylarginine is elevated in asymptomatic, untreated HIV-1 infection and correlates with markers of immune activation and disease progression**  
Ann. Clin. Biochem. 2013; :
- Jawalekar, SL; Karnik, A; Bhutey, A  
**Risk of cardiovascular diseases in diabetes mellitus and serum concentration of asymmetrical dimethylarginine**  
Biochem Res Int 2013; 2013: 189430
- Kleparnik, M; Tomandlova, M; Glatz, Z; Tomandl, J  
**Determination of asymmetric and symmetric dimethylarginines in human plasma by HPLC with electrochemical detection**  
J Sep Sci 2013; 36: 3696-3701
- Korkosz, M; Gaşowski, J; Surdacki, A; Leszczyński, P; Pawlak-Buś, K; Jeka, S; Siedlar, M; Grodzicki, T  
**Disparate effects of anti-TNF- $\alpha$  therapies on measures of disease activity and mediators of endothelial damage in ankylosing spondylitis**  
Pharmacol Rep 2013; 65: 891-897
- Kraml, PJ; Syrovátka, P; Potočková, J; Anděl, M  
**The oxidized low-density lipoprotein/ $\beta$ 2-glycoprotein I complex is associated with abdominal obesity in healthy middle-aged men**  
Ann. Nutr. Metab. 2013; 62:
- Kruszelnicka, O; Surdacki, A; Golay, A  
**Differential associations of angiographic extent and severity of coronary artery disease with asymmetric dimethylarginine but not insulin resistance in non-diabetic men with stable angina: a cross-sectional study**  
Cardiovasc Diabetol 2013; 12: 145
- Labrujere, S; van Houten, EL; de Vries, R; Musterd-Bagghoe, UM; Garrelds, IM; Kramer, P; Danser, AH; Villalón, CM; Visser, JA; Van Den Brink, AM  
**Analysis of the vascular responses in a murine model of polycystic ovary syndrome**  
J. Endocrinol. 2013; 218: 205-213
- Liao, YC; Wang, YS; Guo, YC; Lin, WL; Chang, MH; Hank Juo, SH  
**Let-7g improves multiple endothelial functions through targeting TGF- $\beta$  and SIRT-1 signaling**  
J. Am. Coll. Cardiol. 2013; :
- Miralbell, J; López-Cancio, E; López-Oloriz, J; Arenillas, JF; Barrios, M; Soriano-Raya, JJ; Galán, A; Cáceres, C; Alzamora, M; Pera, G; Toran, P; Dávalos, A; Mataró, M  
**Cognitive patterns in relation to biomarkers of cerebrovascular disease and vascular risk factors**  
Cerebrovasc. Dis. 2013; 36: 98-105
- Nagasaka, H; Miida, T; Yorifuji, T; Hirano, K; Inui, A; Fujisawa, T; Tsukahara, H; Hayashi, H; Inomata, Y  
**Metabolic improvements in intrahepatic porto-systemic venous shunt presenting various metabolic abnormalities by 4-phenylacetate**  
Clin. Chim. Acta 2013; 419: 52-56
- Nagasaka, H; Yorifuji, T; Egawa, H; Inui, A; Fujisawa, T; Komatsu, H; Tsukahara, H; Uemoto, S; Inomata, Y  
**Characteristics of NO cycle coupling with urea cycle in non-hyperammonemic carriers of ornithine transcarbamylase deficiency**  
Mol. Genet. Metab. 2013; 109: 251-264
- Noorbakhsh, M; Kianpour, M; Nematbakhsh, M  
**Serum levels of asymmetric dimethylarginine, vascular endothelial growth factor, and nitric oxide metabolite levels in preeclampsia patients**  
ISRN Obstet Gynecol 2013; 2013: 104213
- Pamuk, BO; Sari, I; Selcuk, S; Gokce, G; Kozaci, DL  
**Evaluation of circulating endothelial biomarkers in familial Mediterranean fever**  
Rheumatol. Int. 2013; 33: 1967-1972
- Ras, RT; Zock, PL; Zebregs, YE; Johnston, NR; Webb, DJ; Draijer, R  
**Effect of polyphenol-rich grape seed extract on ambulatory blood pressure in subjects with pre-and stage I hypertension**  
Br. J. Nutr. 2013; 110: 2234-2241
- Shokouh, P; Joharimoghadam, A; Roohafza, H; Sadeghi, M; Golabchi, A; Boshtam, M; Sarrafzadegan, N  
**Effects of Pioglitazone on Asymmetric Dimethylarginine and Components of the Metabolic Syndrome in Nondiabetic Patients (EPICAMP Study): A Double-Blind, Randomized Clinical Trial**  
PPAR Res 2013; 2013:

- Spasovski, D; Latifi, A; Osmani, B; Krstevska-Balkanov, S; Kafedizska, I; Slaninka-Micevska, M; Dejanova, B; Alabakovska, S; Balkanov, T  
**Determination of the diagnostic values of asymmetric dimethylarginine as an indicator for evaluation of the endothelial dysfunction in patients with rheumatoid arthritis**  
Arthritis 2013; :
- Surdacki, A; Kruszelnicka, O; Rakowski, T; Jaźwińska-Kozuba, A; Dubiel, JS  
**Asymmetric dimethylarginine predicts decline of glucose tolerance in men with stable coronary artery disease: a 4.5-year follow-up study**  
Cardiovasc Diabetol 2013; 12: 64
- Tomada, N; Tomada, I; Botelho, F; Pacheco-Figueiredo, L; Lopes, T; Negrão, R; Pestana, M; Cruz, F  
**Endothelial function in patients with metabolic syndrome and erectile dysfunction: a question of angiotensin imbalance?**  
Andrology 2013; 1: 541-548
- Witasz, A; Rydén, M; Carrero, JJ; Qureshi, AR; Nordfors, L; Näslund, E; Hammarqvist, F; Arefin, S; Kublickiene, K; Stenvinkel, P  
**Elevated circulating levels and tissue expression of pentraxin 3 in uremia: a reflection of endothelial dysfunction**  
PLoS ONE 2013; 8: e63493
- Yamada, S; Saitoh, S; Machii, H  
**Coronary Artery Spasm Related to Thiol Oxidation and Senescence Marker Protein-30 in Aging**  
Antioxidants & Redox Signaling 2013; 19: 1063-1073
- Yusuf, M; Thaha, M; Yogiarto, RM; Aminuddin, M  
**Dimethylarginine Dimethylaminohydrolase 2 Gene Polymorphism and Its Association with Asymmetrical Dimethyl Arginine in Hemodialyzed Patients**  
Open Journal of Nephrology 2013; 3: 75-81
- Chiang, CH; Huang, PH; Chung, FP; Chen, ZY; Leu, HB; Huang, CC; Wu, TC; Chen, JW; Lin, SJ  
**Decreased circulating endothelial progenitor cell levels and function in patients with nonalcoholic fatty liver disease**  
PLoS ONE 2012; 7: e31799
- Enseleit, F; Sudano, I; Périat, D; Winnik, S; Wolfrum, M; Flammer, AJ; Fröhlich, GM; Kaiser, P; Hirt, A; Haile, SR; Krasniqi, N; Matter, CM; Uhlenhut, K; Högger, P; Neidhart, M; Lüscher, TF; Ruschitzka, F; Noll, G  
**Effects of Pycnogenol on endothelial function in patients with stable coronary artery disease: a double-blind, randomized, placebo-controlled, cross-over study**  
Eur. Heart J. 2012; 33: 1589-1597
- Feldman, L; Abu Hamad, R; Efrati, S; Ashker, A  
**Effect of N-Acetylcysteine on Residual Renal Function in Chronic Haemodialysis Patients Treated with High-Flux Synthetic Dialysis Membranes: A Pilot Study**  
ISRN Nephrology 2012; 2013:
- Feldman, L; Shani, M; Sinuani, I; Beberashvili, I; Weissgarten, J  
**N-acetylcysteine may improve residual renal function in hemodialysis patients: a pilot study**  
Hemodial Int 2012; 16: 512-516
- Frieling, H; Leitmeier, V; Haschemi-Nassab, M; Kornhuber, J; Rhein, M; Bleich, S; Hillemacher, T  
**Reduced plasma levels of asymmetric dimethylarginine (ADMA) in patients with alcohol dependence normalize during withdrawal**  
Eur Neuropsychopharmacol 2012; 22: 836-840
- Huang, CY; Yeh, JS; Shih, CM; Lin, FY  
**The Association between Serum Asymmetric Dimethylarginine and Coronary Atherosclerotic Plaque in an Asymptomatic Population**  
Immunology, Endocrine & Metabolic Agents in Medicinal Chemistry 2012; 12: 236-243
- Işıklar, OO; Barutcuoğlu, B; Kabaroglu, C; Mutaf, I; Özmen, D; Bayındır, O; Zoghi, M; Uluer, H  
**Do cardiac risk factors affect the homocysteine and asymmetric dimethylarginine relationship in patients with coronary artery diseases?**  
Clin. Biochem. 2012; 45: 1325-1330
- Ivarsen, P; Povlsen, JV; Christensen, KL  
**Effect of alfalcidol on cardiac function in patients with chronic kidney disease stage 4 and secondary hyperparathyroidism: a pilot study**  
Scand. J. Urol. Nephrol. 2012; 46: 381-388
- Jung, CS; Lange, B; Zimmermann, M; Seifert, V  
**The CSF concentration of ADMA, but not of ET-1, is correlated with the occurrence and severity of cerebral vasospasm after subarachnoid hemorrhage**  
Neurosci. Lett. 2012; 524: 20-24

## 2012

- Alsagaff, MY; Thaha, M; Aminuddin, M; Yogiarto, RM; Yogiartoro, M; Tomino, Y  
**Asymmetric dimethylarginine: a novel cardiovascular risk factor in end-stage renal disease**  
J. Int. Med. Res. 2012; 40: 340-349
- Aydin, M; Koca, C; Uysal, S; TOTAN, Y  
**Serum nitric oxide, asymmetric dimethylarginine, and plasma homocysteine levels in active Behçet's disease**  
Turk J Med Sci 2012; 42: 1194-1199
- Briceño, LG; Alfieria, AB; Briceño, AB  
**Los niveles plasmáticos de citocinas inflamatorias están elevados en pacientes con insuficiencia cardiaca**  
Revista Latinoamericana de Hipertension 2012; 4:

- Kadrabova, J; Krajcovicova-Kudlackova, M  
**Protective properties of complex of quercetin, selenium, catechins and curcumin against DNA damage**  
Oxidants and Antioxidants in Medical Science 2012; 1: 179
- Khan, BV; Rahman, ST; Haque, T; Merchant, N; Bhaheetharan, S; Harris, J; Umar, K; Wahi, J; Ferdinand, KC;  
**Vascular effects of nebivolol added to hydrochlorothiazide in African Americans with hypertension and echocardiographic evidence of diastolic dysfunction: the NASAA study**  
J. Cardiovasc. Pharmacol. Ther. 2012; 17: 291-307
- Khalifa, NM; Gad, MZ; Hataba, AA; Mahran, LG  
**Changes in ADMA and TAFI levels after stenting in coronary artery disease patients**  
Mol Med Rep 2012; 6: 855-859
- Khazaei, M; Zarei, M; Esmaeil, N  
**Serum Asymmetric Dimethylarginine Concentrations in Two Models of Experimentally Induced Hypertension**  
ISRN Nephrology 2012; :
- King, DE; Player, M; Everett, CJ  
**The impact of pioglitazone on ADMA and oxidative stress markers in patients with type 2 diabetes**  
Prim Care Diabetes 2012; 6: 157-161
- Korkosz, M; Sulicka, J; Surdacki, A  
**Dimethylated L-arginine analogues versus autoantibodies in early rheumatoid arthritis**  
Scand. J. Rheumatol. 2012; 41: 82-83
- Li, J; Xia, W; Feng, W; Qu, X  
**Effects of rosuvastatin on serum asymmetric dimethylarginine levels and atrial structural remodeling in atrial fibrillation dogs**  
Pacing Clin Electrophysiol 2012; 35: 456-464
- López-Cancio, E; Galán, A; Dorado, L; Jiménez, M; Hernández, M; Millán, M; Reverté, S; Suñol, A; Barallat, J; Massuet, A; Alzamora, MT; Dávalos, A; Arenillas, JF  
**Biological signatures of asymptomatic extra- and intracranial atherosclerosis: the Barcelona-AsIA (Asymptomatic Intracranial Atherosclerosis) study**  
Stroke 2012; 43: 2712-2719
- Luksha, L; Stenvinkel, P; Hammarqvist, F; Carrero, JJ; Davidge, ST; Kublickiene, K  
**Mechanisms of endothelial dysfunction in resistance arteries from patients with end-stage renal disease**  
PLoS ONE 2012; 7: e36056
- Moreno-Luna, R; Muñoz-Hernandez, R; Miranda, ML; Costa, AF; Jimenez-Jimenez, L; Vallejo-Vaz, AJ; Muriana, FJ; Villar, J; Stiefel, P  
**Olive oil polyphenols decrease blood pressure and improve endothelial function in young women with mild hypertension**  
Am. J. Hypertens. 2012; 25: 1299-1304
- Nakhjavani, M; Morteza, A; Asgarani, F; Khalilzadeh, O; Ghazizadeh, Z; Bathaie, SZ; Esteghamati, A  
**The dual behavior of heat shock protein 70 and asymmetric dimethylarginine in relation to serum CRP levels in type 2 diabetes**  
Gene 2012; 498: 107-111
- Napora, M; Graczykowska, A; Próchniewska, K; Zdrojewski, Z; Całka, A; Górny, J; Stompór, T  
**Relationship between serum asymmetric dimethylarginine and left ventricular structure and function in patients with end-stage renal disease treated with hemodialysis**  
Pol. Arch. Med. Wewn. 2012; 122: 226-234
- Parenica, J; Nemeč, P; Tomandl, J; Ondrasek, J; Pavkova-Goldbergova, M; Tretina, M; Jarkovsky, J; Littnerova, S; Poloczek, M; Pokorný, P; Spinar, J; Cermakova, Z; Miklik, R; Malík, P; Pes, O; Lipkova, J; Tomandlova, M; Kala, P  
**Prognostic utility of biomarkers in predicting of one-year outcomes in patients with aortic stenosis treated with transcatheter or surgical aortic valve implantation**  
PLoS ONE 2012; 7: e48851
- Rueda-Clausen, CF; Córdoba-Porras, A; Bedoya, G; Silva, FA; Zarruk, JG; López-Jaramillo, P; Villa, LA  
**Increased plasma levels of total homocysteine but not asymmetric dimethylarginine in Hispanic subjects with ischemic stroke FREC-VI sub-study**  
Eur. J. Neurol. 2012; 19: 417-425
- Salum, E; Kampus, P; Zilmer, M; Eha, J; Butlin, M; Avolio, AP; Põdräma, T; Arend, A; Aunapuu, M; Kals, J  
**Effect of vitamin D on aortic remodeling in streptozotocin-induced diabetes**  
Cardiovasc Diabetol 2012; 11: 58
- Serg, M; Kampus, P; Kals, J; Zagura, M; Zilmer, M; Zilmer, K; Kullisaar, T; Eha, J  
**Nebivolol and metoprolol: long-term effects on inflammation and oxidative stress in essential hypertension**  
Scand. J. Clin. Lab. Invest. 2012; 72: 427-432
- Sulicka, J; Surdacki, A; Strach, M; Kwater, A; Gryglewska, B; Œwiklińska, M; Balwierz, W; Grodzicki, TK  
**Elevated asymmetric dimethylarginine in young adult survivors of childhood acute lymphoblastic leukemia: a preliminary report**  
Dis. Markers 2012; 33: 69-76
- Sztanek, F; Seres, I; Harangi, M; Lőcsey, L; Koncsos, P; Paragh, G  
**Effect of nutritional status on human paraoxonase-1 activity in patients with chronic kidney disease**  
Kidney Blood Press. Res. 2012; 36: 310-319

- Sztanek, F; Seres, I; Harangi, M; Lőcsey, L; Padra, J; Paragh, GJ; Asztalos, L; Paragh, G  
**Decreased paraoxonase 1 (PON1) lactonase activity in hemodialyzed and renal transplanted patients. A novel cardiovascular biomarker in end-stage renal disease**  
Nephrol. Dial. Transplant. 2012; 27: 2866-2872
- Teplan, V; Racek, J  
**Asymmetric Dimethylarginine, in Uremic Toxins (ed T. Niwa)**  
Uremic Toxins 2012; :
- Thomson, RL; Brinkworth, GD; Noakes, M; Clifton, PM; Norman, RJ; Buckley, JD  
**The effect of diet and exercise on markers of endothelial function in overweight and obese women with polycystic ovary syndrome**  
Hum. Reprod. 2012; 27: 2169-2176
- Vladimirova-Kitova, LG; Deneva-Koycheva, TI  
**The effect of simvastatin on asymmetric dimethylarginine and flow-mediated vasodilation after optimizing the LDL level: a randomized, placebo-controlled study**  
Vascul. Pharmacol. 2012; 56: 122-130
- Wibowo, N; Purwosunu, Y; Sekizawa, A; Farina, A; Idriansyah, L; Fitriana, I  
**Antioxidant supplementation in pregnant women with low antioxidant status**  
J. Obstet. Gynaecol. Res. 2012; 38: 1152-1161
- Eloit, S; Schepers, E; Barreto, DV; Barreto, FC; Liabeuf, S; Van Biesen, W; Verbeke, F; Glorieux, G; Choukroun, G; Massy, Z; Vanholder, R  
**Estimated glomerular filtration rate is a poor predictor of concentration for a broad range of uremic toxins**  
Clin J Am Soc Nephrol 2011; 6: 1266-1273
- Erre, GL; Sanna, P; Zinellu, A; Ponchietti, A; Fenu, P; Sotgia, S; Carru, C; Ganau, A; Passiu, G  
**Plasma asymmetric dimethylarginine (ADMA) levels and atherosclerotic disease in ankylosing spondylitis: a cross-sectional study**  
Clin. Rheumatol. 2011; 30: 21-7
- Grassi, G; Seravalle, G; Ghiadoni, L; Tripepi, G; Bruno, RM; Mancina, G; Zoccali, C  
**Sympathetic nerve traffic and asymmetric dimethylarginine in chronic kidney disease**  
Clin J Am Soc Nephrol 2011; 6: 2620-2627
- Haghikia, A; Missol-Kolka, E; Tsikas, D; Venturini, L; Brundiers, S; Castoldi, M; Muckenthaler, MU; Eder, M; Stapel, B; Thum, T; Haghikia, A; Petrasch-Parwez, E; Drexler, H; Hilfiker-Kleiner, D; Scherr, M  
**Signal transducer and activator of transcription 3-mediated regulation of miR-199a-5p links cardiomyocyte and endothelial cell function in the heart: a key role for ubiquitin-conjugating enzymes**  
Eur. Heart J. 2011; 32: 1287-1297
- Hanssen, H; Nickel, T; Drexler, V; Hertel, G; Emslander, I; Sisic, Z; Lorang, D; Schuster, T; Kotliar, KE; Pressler, A; Schmidt-Trucksäss, A; Weis, M; Halle, M  
**Exercise-induced alterations of retinal vessel diameters and cardiovascular risk reduction in obesity**  
Atherosclerosis 2011; 216: 433-439
- Hermansen, SE; Kalstad, T; How, OJ; Myrnes, T  
**Inflammation and reduced endothelial function in the course of severe acute heart failure**  
Transl Res 2011; 157: 117-127
- Hu, X; Atzler, D; Xu, X; Zhang, P; Guo, H; Lu, Z; Fassett, J; Schwedhelm, E; Böger, RH; Bache, RJ; Chen, Y  
**Dimethylarginine dimethylaminohydrolase-1 is the critical enzyme for degrading the cardiovascular risk factor asymmetrical dimethylarginine**  
Arterioscler. Thromb. Vasc. Biol. 2011; 31: 1540-1546
- Ioakeimidis, N; Vlachopoulos, C; Rokkas, K; Aggelis, A; Terentes-Printzios, D; Samentzas, A; Alexopoulos, N; Stefanadis, C  
**Relationship of asymmetric dimethylarginine with penile Doppler ultrasound parameters in men with vasculogenic erectile dysfunction**  
Eur. Urol. 2011; 59: 948-955

## 2011

- Andrade, F; Rodríguez-Soriano, J; Prieto, JA; Aguirre, M; Ariceta, G; Lage, S; Azcona, I; Prado, C; Sanjurjo, P; Aldámiz-Echevarría, L  
**Methylation cycle, arginine-creatinine pathway and asymmetric dimethylarginine in paediatric renal transplant**  
Nephrol. Dial. Transplant. 2011; 26: 328-236
- Carlsson, M; Lindström, T; Nyström, F  
**Evidence of Daytime Variation of Asymmetric Dimethylarginine: An Intervention Study with Rosiglitazone in Type 2 Diabetes**  
The Open Endocrinology Journal 2011; 5: 14-18
- Damiati, SA; Khoja, SM  
**Serum Asymmetric Dimethyl-L-Arginine in Renal Failure Patients Living in Jeddah Region, Saudi Arabia**  
Trends in Medical Research 2011; 6: 14-22
- Deneva-Koycheva, TI; Vladimirova-Kitova, LG; Angelova, EA; Tsvetkova, TZ  
**Plasma asymmetric dimethylarginine levels in healthy people**  
Folia Med (Plovdiv) 2011; 53: 28-33

- Kilic, S; Yilmaz, N; Zulfikaroglu, E; Erdogan, G; Aydin, M; Batioglu, S  
**Inflammatory-metabolic parameters in obese and nonobese normoandrogenemic polycystic ovary syndrome during metformin and oral contraceptive treatment**  
Gynecol. Endocrinol. 2011; 27: 622-629
- Kocak, H; Oner-Iyidogan, Y; Gurdol, F; Oner, P; Esin, D  
**Serum asymmetric dimethylarginine and nitric oxide levels in obese postmenopausal women**  
J. Clin. Lab. Anal. 2011; 25: 174-178
- Koncsos, P; Seres, I; Harangi, M; Páll, D; Józsa, L; Bajnok, L; Nagy, EV; Paragh, G  
**Favorable effect of short-term lifestyle intervention on human paraoxonase-1 activity and adipokine levels in childhood obesity**  
J Am Coll Nutr 2011; 30: 333-339
- Kuncová, J; Chvojka, J; Sýkora, R; Svíglerová, J; Stengl, M; Nalos, L; Kroužecký, A; Matějovič, M  
**Tissue concentrations of vasoactive intestinal peptide are affected by peritonitis-induced sepsis and hemofiltration in pigs**  
Physiol Res 2011; 60: 531-540
- Kuo, KL; Hung, SC; Tarng, DC; Selim, G  
**Relationship between endothelial adhesion of peripheral blood mononuclear cells (pbmcs) and superoxide production and of inflammatory markers in hemodialysis (hd) patients**  
Ndt plus 2011; 4:
- La Favor, JD; Hollis, BC; Mokshagundam, SL; Olive, JL  
**Serum hsCRP and visfatin are elevated and correlate to carotid arterial stiffness in spinal cord-injured subjects**  
Spinal Cord 2011; 49: 961-966
- Li Volti, G; Salomone, S; Sorrenti, V; Mangiameli, A; Urso, V; Siarkos, I; Galvano, F; Salamone, F  
**Effect of silibinin on endothelial dysfunction and ADMA levels in obese diabetic mice**  
Cardiovasc Diabetol 2011; 10: 62
- Lüersen, K; Schmelzer, C; Boesch-Saadatmandi, C; Kohl, C; Rimbach, G; Döring, F  
**Paraoxonase 1 polymorphism Q192R affects the pro-inflammatory cytokine TNF-alpha in healthy males**  
BMC Res Notes 2011; 4: 141
- Luksha, N; Luksha, L; Carrero, JJ; Hammarqvist, F; Stenvinkel, P; Kublickiene, K  
**Impaired resistance artery function in patients with end-stage renal disease**  
Clin. Sci. 2011; 120: 525-536
- Mihout, F; Shweke, N; Bigé, N; Jouanneau, C; Dussault, JC; Ronco, P; Chatziantoniou, C; Boffa, JJ  
**Asymmetric dimethylarginine (ADMA) induces chronic kidney disease through a mechanism involving collagen and TGF-β1 synthesis**  
J. Pathol. 2011; 223: 37-45
- Puls, M; Schroeter, MR; Steier, J; Stijohann, L; Hasenfuss, G; Konstantinides, S; Schäfer, K  
**Effect of smoking cessation on the number and adhesive properties of early outgrowth endothelial progenitor cells**  
Int. J. Cardiol. 2011; 152: 61-69
- Rossenbach, J; Mueller, GA; Lange, K; Armstrong, VW; Schmitto, JD; Hintze, E; Helfmann, J; Konstantinides, S; Koziolok, MJ  
**Lipid-apheresis improves microcirculation of the upper limbs**  
J Clin Apher 2011; 26: 167-173
- Sanayama, Y; Nagasaka, H; Takayanagi, M; Ohura, T; Sakamoto, O; Ito, T; Ishige-Wada, M; Usui, H; Yoshino, M; Ohtake, A; Yorifuji, T; Tsukahara, H; Hirayama, S; Miida, T; Fukui, M; Okano, Y  
**Experimental evidence that phenylalanine is strongly associated to oxidative stress in adolescents and adults with phenylketonuria**  
Mol. Genet. Metab. 2011; 103: 220-225
- Schepers, E; Barreto, DV; Liabeuf, S; Glorieux, G; Eloit, S; Barreto, FC; Massy, Z; Vanholder, R; European Uremic Toxin Work Group (EUTox),  
**Symmetric dimethylarginine as a proinflammatory agent in chronic kidney disease**  
Clin J Am Soc Nephrol 2011; 6: 2374-2383
- Schlager, O; Giurgea, A; Schuhfried, O; Seidinger, D; Hammer, A; Gröger, M; Fialka-Moser, V; Gschwandtner, M; Koppensteiner, R; Steiner, S  
**Exercise training increases endothelial progenitor cells and decreases asymmetric dimethylarginine in peripheral arterial disease: a randomized controlled trial**  
Atherosclerosis 2011; 217: 240-248
- Serg, M; Kampus, P; Kals, J; Zagura, M; Muda, P; Tuomainen, TP; Zilmer, K; Salum, E; Zilmer, M; Eha, J  
**Association between asymmetric dimethylarginine and indices of vascular function in patients with essential hypertension**  
Blood Press. 2011; 20: 111-116
- Sun, L; Zhang, T; Yu, X; Xin, W; Lan, X; Zhang, D; Huang, C; Du, G  
**Asymmetric dimethylarginine confers the communication between endothelial and smooth muscle cells and leads to VSMC migration through p38 and ERK1/2 signaling cascade**  
FEBS Lett. 2011; 585: 2727-2734
- Topal, G; Donmez, A; Doğan, BS; Kucur, M; Cengiz, DT; Berkoz, FB; Erdogan, N  
**Asymmetric dimethylarginine (ADMA) levels are increased in patients with fibromyalgia: correlation with tumor necrosis factor-α (TNF-α) and 8-iso-prostaglandin F(2α) (8-iso-PGF(2α))**  
Clin. Biochem. 2011; 44: 364-367



- Tousoulis, D; Bouras, G; Antoniadis, C; Marinou, K; Papageorgiou, N; Miliou, A; Hatzis, G; Stefanadi, E; Tsioufis, C; Stefanadis, C  
**Methionine-induced homocysteinemia impairs endothelial function in hypertensives: the role of asymmetrical dimethylarginine and antioxidant vitamins**  
Am. J. Hypertens. 2011; 24: 936-942
- Tuygun, AK; Tuygun, A; Yurtseven, N; Şensöz, Y  
**Taux de diméthylarginine asymétrique au cours de la maladie de Buerger**  
Annales de Chirurgie Vasculaire 2011; 25: 589-596
- Tuygun, AK; Tuygun, A; Yurtseven, N; Sensöz, Y; Günay, R; Keser, M; Tuygun, UH; Ketenci, B; Sahin, S; Yekeler, I  
**Asymmetric dimethylarginine levels in Buerger's disease**  
Ann Vasc Surg 2011; 25: 547-554
- Vladimirova-Kitova, LG; Deneva, TI; Marinov, B  
**Effect of moderate and high-dose simvastatin on asymmetric dimethylarginine-homocysteine metabolic pathways in patients with newly detected severe hypercholesterolemia**  
Cardiovasc Ther 2011; 29: 340-348
- Vladimirova-Kitova, LG; Deneva-Koicheva, TI  
**Increased intima-media thickness in carriers of the ldl-receptor defective gene versus noncarriers with newly detected asymptomatic severe hypercholesterolemia**  
Echocardiography 2011; 28: 223-234
- Wang, Y; Zhang, M; Liu, Y; Liu, Y; Chen, M  
**The effect of nebivolol on asymmetric dimethylarginine system in spontaneously hypertension rats**  
Vascul. Pharmacol. 2011; 54: 36-43
- Böger, RH; Hilgers, KF; Jacobi, J  
**Role of asymmetric dimethylarginine for angiotensin for angiotensin II-induced target organ damage in mice**  
Am J Physiol Heart Circ Physiol 2010; 294: 1058-1066
- Cavusoglu, E; Ruwende, C; Chopra, V; Poludasu, S; Yanamadala, S; Frishman, WH; Eng, C; Pinsky, DJ; Marmur, JD  
**Relation of baseline plasma ADMA levels to cardiovascular morbidity and mortality at two years in men with diabetes mellitus referred for coronary angiography**  
Atherosclerosis 2010; 210: 226-231
- Dimitroulas, T; Giannakoulas, G; Papadopoulou, K; Karvounis, H; Dimitroula, H; Koliakos, G; Karamitsos, T; Parcharidou, D; Settas, L  
**Early detection of cardiac involvement in systemic sclerosis assessed by tissue-Doppler echocardiography: relationship with neurohormonal activation and endothelial dysfunction**  
J. Rheumatol. 2010; 37: 993-999
- Dimitroulas, T; Giannakoulas, G; Papadopoulou, K; Sfetsios, T; Karvounis, H; Dimitroula, H; Parcharidou, D; Koliakos, G; Garyfallos, A; Styliadis, I; Settas, L  
**Left atrial volume and N-terminal pro-B type natriuretic peptide are associated with elevated pulmonary artery pressure in patients with systemic sclerosis**  
Clin. Rheumatol. 2010; 29: 957-964
- Eiselt, J; Rajdl, D; Racek, J; Siroká, R; Trefil, L; Opatrná, S  
**Asymmetric dimethylarginine in hemodialysis, hemodiafiltration, and peritoneal dialysis**  
Artif Organs 2010; 34: 420-425
- Galyavich, A; Khamidullina, A; Galyavi, R  
**Effects of Antihypertensive Drugs on Plasma Levels of Nitric Oxide Derivates Products and Asymmetric Dimethylarginine in Target Achieved Systolic Blood Pressure Patients: Pp.6.263**  
Journal of Hypertension 2010; 28: 128

## 2010

- Ames, PRJ; Ortiz-Cadenas, A  
**Rosuvastatin treatment is associated with a decrease of serum oxidised low-density lipoprotein/beta2-glycoprotein I complex concentration in type 2 diabetes**  
The British Journal of Diabetes & Vascular Disease 2010; 10: 292-299
- An, H; An, S; Erdoğan, E; Tiryakioğlu, O; Huysal, K; Koca, V; Bozat, T  
**The effects of endothelial dysfunction and inflammation on slow coronary flow**  
Turk Kardiyol Dern Ars 2010; 38: 327-333
- Ari, H; Ari, S; Erdoğan, E; Tiryakioğlu, O; Ustündağ, Y; Huysal, K; Koca, V; Bozat, T  
**A novel predictor of restenosis and adverse cardiac events: asymmetric dimethylarginine**  
Heart Vessels 2010; 25: 19-26
- Głowińska-Olszewska, B; Luczyński, W; Jabłońska, J; Otocka, A; Florys, B; Bossowski, A  
**Asymmetric dimethylarginine (ADMA) in children with diabetes type 1**  
Pediatri Endocrinol Diabetes Metab 2010; 16: 137-141
- Gurdol, F; Cimsit, M; Oner-Iyidogan, Y; Kocak, H; Sengun, S; Yalcinkaya-Demirsoz, S  
**Collagen synthesis, nitric oxide and asymmetric dimethylarginine in diabetic subjects undergoing hyperbaric oxygen therapy**  
Physiol Res 2010; 59: 423-429

- Heida, NM; Müller, JP; Cheng, IF; Leifheit-Nestler, M; Faustin, V; Riggert, J; Hasenfuss, G; Konstantinides, S; Schäfer, K  
**Effects of obesity and weight loss on the functional properties of early outgrowth endothelial progenitor cells**  
J. Am. Coll. Cardiol. 2010; 55: 357-367
- Kilic, S; Yilmaz, N; Erdogan, G; Aydin, M; Tasdemir, N; Doganay, M; Batioglu, S  
**Effect of non-oral estrogen on risk markers for metabolic syndrome in early surgically menopausal women**  
Climacteric 2010; 13: 55-62
- Kimoto, M; Bache, RJ; Chen, YJ; Li, Y; Zhang, P  
**Dimethylarginine dimethylaminohydrolase and endothelial dysfunction in failing hearts**  
Am J Physiol Heart Circ Physiol 2010; 289: 2212-2219
- Lee, KH; Gil, HW; Yang, JO; Lee, EY; Hong, SY  
**Sequential change of asymmetric dimethylarginine levels after initiation of peritoneal dialysis in patients with end-stage renal disease**  
Nephrology (Carlton) 2010; 15: 639-643
- Meng, J; Li, ZX; Jiang, W; Xu, C; Li, YC  
**Relationship of Serum ADMA With pulmonary hypertension in patients on hemodialysis**  
Dialysis & Transplantation 2010; 39: 242-246
- Mohamadin, AM; Habib, FA; Al-Saggaf, AA  
**Cardiovascular disease markers in women with polycystic ovary syndrome with emphasis on asymmetric dimethylarginine and homocysteine**  
Ann Saudi Med 2010; 30: 278-283
- Nagasaka, H; Okano, Y; Aizawa, M; Miida, T; Yorifuji, T; Tajima, G; Sakura, N; Takatani, T; Sanayama, Y; Sugamoto, K; Mayumi, M; Kobayashi, K; Hirano, K; Takayanagi, M; Tsukahara, H  
**Altered metabolisms of mediators controlling vascular function and enhanced oxidative stress in asymptomatic children with congenital portosystemic venous shunt**  
Metab. Clin. Exp. 2010; 59: 107-113
- Nakhjavani, M; Karimi-Jafari, H; Esteghamati, A; Khalilzadeh, O; Asgarani, F; Ghadiri-Anari, A  
**ADMA is a correlate of insulin resistance in early-stage diabetes independent of hs-CRP and body adiposity**  
Ann. Endocrinol. (Paris) 2010; 71: 303-308
- Napoli, R; Guardasole, V; Zarra, E; D'Anna, C  
**Impaired endothelial- and nonendothelial-mediated vasodilation in patients with acute or chronic hypothyroidism**  
Clinical Endocrinology 2010; 72: 107-111
- Nascimento, MM; Suliman, ME; Silva, M; Chinaglia, T; Marchioro, J; Hayashi, SY; Riella, MC; Lindholm, B; Anderstam, B  
**Effect of oral N-acetylcysteine treatment on plasma inflammatory and oxidative stress markers in peritoneal dialysis patients: a placebo-controlled study**  
Perit Dial Int 2010; 30: 336-342
- Perna, M; Roman, MJ; Alpert, DR; Crow, MK; Lockshin, MD; Sammaritano, L; Devereux, RB; Cooke, JP; Salmon, JE  
**Relationship of asymmetric dimethylarginine and homocysteine to vascular aging in systemic lupus erythematosus patients**  
Arthritis Rheum. 2010; 62: 1718-1722
- Quinaglia, T; Figueiredo, VN; Martins, LC  
**Endothelial Dysfunction in Refractory Hypertension: Correlation Between Blood Pressure Levels and Nocturnal Dipping Bp Impairment: Pp. 6.261**  
Journal of Hypertension 2010; 28: 127-128
- Sandrim, VC; Palei, AC; Metzger, IF; Cavalli, RC; Duarte, G; Tanus-Santos, JE  
**Interethnic differences in ADMA concentrations and negative association with nitric oxide formation in preeclampsia**  
Clin. Chim. Acta 2010; 411: 1457-1460
- Santilli, F; Formoso, G; Sbraccia, P; Averna, M; Miccoli, R; Di Fulvio, P; Ganci, A; Pulizzi, N; Lattanzio, S; Ciabattone, G; Consoli, A; Lauro, R; Patrono, C; Davì, G  
**Postprandial hyperglycemia is a determinant of platelet activation in early type 2 diabetes mellitus**  
J. Thromb. Haemost. 2010; 8: 828-837
- Savoia, C; Ebrahimian, T; Lemarié, CA; Paradis, P; Iglarz, M; Amiri, F; Javeshgani, D; Schiffrin, EL  
**Countervailing vascular effects of rosiglitazone in high cardiovascular risk mice: role of oxidative stress and PRMT-1**  
Clin. Sci. 2010; 118: 583-592
- Shakoor, SK; Aldibbiat, A; Ingoe, LE; Campbell, SC; Sibal, L; Shaw, J; Home, PD; Razvi, S; Weaver, JU  
**Endothelial progenitor cells in subclinical hypothyroidism: the effect of thyroid hormone replacement therapy**  
J. Clin. Endocrinol. Metab. 2010; 95: 319-322
- Surdacki, A; Marewicz, E; Wieczorek-Surdacka, E; Rakowski, T; Szastak, G; Pryjma, J; Dudek, D; Dubiel, JS  
**Synergistic effects of asymmetrical dimethyl-L-arginine accumulation and endothelial progenitor cell deficiency on renal function decline during a 2-year follow-up in stable angina**  
Nephrol. Dial. Transplant. 2010; 25: 2576-2583

- Sydow, K; Fortmann, SP; Fair, JM; Varady, A; Hlatky, MA; Go, AS; Iribarren, C; Tsao, PS; ADVANCE Investigators,

**Distribution of asymmetric dimethylarginine among 980 healthy, older adults of different ethnicities**

Clin. Chem. 2010; 56: 111-120

- Teede, HJ; Meyer, C; Hutchison, SK; Zoungas, S; McGrath, BP; Moran, LJ  
**Endothelial function and insulin resistance in polycystic ovary syndrome: the effects of medical therapy**  
Fertil. Steril. 2010; 93: 184-191
- Teplan, V; Vyhnánek, F; Gürlich, R; Haluzík, M; Racek, J; Vyhnankova, I; Stollová, M; Teplan, V  
**Increased proinflammatory cytokine production in adipose tissue of obese patients with chronic kidney disease**  
Wien. Klin. Wochenschr. 2010; 122: 466-473
- Tsioufis, C; Dimitriadis, K; Andrikou, E; Thomopoulos, C; Tsiachris, D; Stefanadi, E; Mihas, C; Miliou, A; Papademetriou, V; Stefanadis, C  
**ADMA, C-reactive protein, and albuminuria in untreated essential hypertension: a cross-sectional study**  
Am. J. Kidney Dis. 2010; 55: 1050-1059
- Tsioufis, C; Thomopoulos, C; Soldatos, N; Syrseloudis, D; Kasiakogias, A; Silvestros, S; Stefanadi, E; Mostratou, E; Stefanadis, C  
**The conjoint detrimental effect of chronic periodontal disease and systemic inflammation on asymmetric dimethyl-arginine in untreated hypertensive subjects**  
Atherosclerosis 2010; 208: 258-263
- Vladimirova-Kitova, L; Deneva, T; Marinov, B  
**Predictors of the intima-media thickness of carotid artery in asymptomatic newly detected severe hypercholesterolemic patients**  
Clin Physiol Funct Imaging 2010; 30: 250-259
- Vladimirova-Kitova, LG; Deneva-Koycheva, T  
**Asymmetric dimethylarginine--a determinant of the effect of the high dose Simvastatin**  
Clin. Biochem. 2010; 43: 843-850
- Vladimirova-Kitova, LG; Terzieva, DD  
**Asymmetric dimethylarginine determines the effect of simvastatin on endothelium-dependent vasodilation in severe hypercholesterolemia**  
Clin.Lipidol 2010; 5: 29-41

## 2009

---

- Antoniadou, C; Shirodaria, C; Leeson, P; Antonopoulos, A; Warrick, N; Van-Assche, T; Cunnington, C; Tousoulis, D; Pillai, R; Ratnatunga, C; Stefanadis, C; Channon, KM  
**Association of plasma asymmetrical dimethylarginine (ADMA) with elevated vascular superoxide production and endothelial nitric oxide synthase uncoupling: implications for endothelial function in human atherosclerosis**  
Eur. Heart J. 2009; 30: 1142-1150
- Assal, HS; Younes, K; Alsayed, A; Hasan, N; Ali, AR  
**Asymmetric Dimethylarginine (ADMA) and Progression of Nephropathy in Patients with Type 2 Diabetes**  
Kidney 2009; 18: 247-251
- Böger, RH; Maas, R; Schulze, F; Schwedhelm, E  
**Asymmetric dimethylarginine (ADMA) as a prospective marker of cardiovascular disease and mortality--an update on patient populations with a wide range of cardiovascular risk**  
Pharmacol. Res. 2009; 60: 481-487
- Cavusoglu, E; Ruwende, C; Chopra, V; Yanamadala, S; Eng, C; Pinsky, DJ; Marmur, JD  
**Relationship of baseline plasma ADMA levels to cardiovascular outcomes at 2 years in men with acute coronary syndrome referred for coronary angiography**  
Coron. Artery Dis. 2009; 20: 112-117
- Chen, Y; Xu, X; Sheng, M; Zhang, X; Gu, Q; Zheng, Z  
**PRMT-1 and DDAHs-induced ADMA upregulation is involved in ROS- and RAS-mediated diabetic retinopathy**  
Exp. Eye Res. 2009; 89: 1028-1034
- Coen, G; Mantella, D; Sardella, D; Beraldi, MP; Ferrari, I; Pierantozzi, A; Lippi, B; Di Giulio, S  
**Asymmetric dimethylarginine, vascular calcifications and parathyroid hormone serum levels in hemodialysis patients**  
J. Nephrol. 2009; 22: 616-622
- Damiati, SA; Khoja, SM  
**Serum concentrations of asymmetric dimethyl-L-arginine in diabetes mellitus and hyperlipidemia**  
Clinical and Experimental Medical Journal 2009; 3: 443-452
- Deshmukh, SH; Patel, SR; Pinassi, E; Mindrescu, C; Hermance, EV; Infantino, MN; Coppola, JT; Staniloae, CS  
**Ranolazine improves endothelial function in patients with stable coronary artery disease**  
Coron. Artery Dis. 2009; 20: 343-347
- Esposito, C; Grosjean, F; Torreggiani, M; Maggi, N; Esposito, V; Migotto, C; Mangione, F; Tinelli, C; Dal Canton, A  
**Increased asymmetric dimethylarginine serum levels are associated with acute rejection in kidney transplant recipients**  
Transplant. Proc. 2009; 41: 1570-1573

- Frank, J; George, TW; Lodge, JK; Rodriguez-Mateos, AM; Spencer, JP; Minihane, AM; Rimbach, G  
**Daily consumption of an aqueous green tea extract supplement does not impair liver function or alter cardiovascular disease risk biomarkers in healthy men**  
J. Nutr. 2009; 139: 58-62
- Guagnano, MT; Ferroni, P; Santilli, F; Paoletti, V; Manigrasso, MR; Pescara, L; Cuccurullo, C; Ciabattoni, G; Davì, G  
**Determinants of platelet activation in hypertensives with microalbuminuria**  
Free Radic. Biol. Med. 2009; 46: 922-927
- Gulhan, I; Kebapcilar, L; Alacacioglu, A; Bilgili, S; Kume, T; Aytac, B; Gunaydin, R  
**Postmenopausal women with osteoporosis may be associated with high endothelin-1**  
Gynecol. Endocrinol. 2009; 25: 674-678
- Hu, X; Xu, X; Zhu, G; Atzler, D; Kimoto, M; Chen, J; Schwedhelm, E; Lüneburg, N; Böger, RH; Zhang, P; Chen, Y  
**Vascular endothelial-specific dimethylarginine dimethylaminohydrolase-1-deficient mice reveal that vascular endothelium plays an important role in removing asymmetric dimethylarginine**  
Circulation 2009; 120: 2222-2229
- Jehlicka, P; Stozický, F; Mayer, O; Varvarovská, J; Racek, J; Trefil, L; Siala, K  
**Asymmetric dimethylarginine and the effect of folate substitution in children with familial hypercholesterolemia and diabetes mellitus type 1**  
Physiol Res 2009; 58: 179-184
- Kebapcilar, L; Yuksel, A; Bozkaya, G; Taner, CE  
**Effects of an EE/CA compared with EE/CA-metformin on serum ADMA levels in women with polycystic ovary syndrome**  
Central European Journal of Medicine 2009; 4: 423-427
- Kuan, Y; Surman, J; Frystyk, J; El Nahas, AM; Flyvbjerg, A; Haylor, JL  
**Lack of effect of IGF-I on the glomerular filtration rate in non-diabetic patients with advanced chronic kidney disease**  
Growth Horm. IGF Res. 2009; 19: 219-225
- Kuncová, J; Svíglerová, J; Kummer, W; Rajdl, D; Chottová-Dvoráková, M; Tonar, Z; Nalos, L; Stengl, M  
**Parasympathetic regulation of heart rate in rats after 5/6 nephrectomy is impaired despite functionally intact cardiac vagal innervation**  
Nephrol. Dial. Transplant. 2009; 24: 2362-2370
- Moran, LJ; Hutchison, SK; Meyer, C; Zoungas, S; Teede, HJ  
**A comprehensive assessment of endothelial function in overweight women with and without polycystic ovary syndrome**  
Clin. Sci. 2009; 116: 761-770
- Nagasaka, H; Okano, Y; Tsukahara, H; Shigematsu, Y; Momoi, T; Yorifuji, J; Miida, T; Ohura, T; Kobayashi, K; Saheki, T; Hirano, K; Takayanagi, M; Yorifuji, T  
**Sustaining hypercitrullinemia, hypercholesterolemia and augmented oxidative stress in Japanese children with aspartate/ glutamate carrier isoform 2-citrin-deficiency even during the silent period**  
Mol. Genet. Metab. 2009; 97: 21-26
- Nagasaka, H; Tsukahara, H; Yorifuji, T; Miida, T; Murayama, K; Tsuruoka, T; Takatani, T; Kanazawa, M; Kobayashi, K; Okano, Y; Takayanagi, M  
**Evaluation of endogenous nitric oxide synthesis in congenital urea cycle enzyme defects.**  
Metabolism Clin. Exp. 2009; 58: 278-282
- Pitocco, D; Zaccardi, F; Di Stasio, E; Romitelli, F; Martini, F; Scaglione, GL; Speranza, D; Santini, S; Zuppi, C; Ghirlanda, G  
**Role of asymmetric-dimethyl-L-arginine (ADMA) and nitrite/nitrate (NOx) in the pathogenesis of oxidative stress in female subjects with uncomplicated type 1 diabetes mellitus**  
Diabetes Res. Clin. Pract. 2009; 86: 173-176
- Puchau, B; Hermsdorff, HH; Zulet, MA; Martínez, JA  
**DDAH2 mRNA expression is inversely associated with some cardiovascular risk-related features in healthy young adults**  
Dis. Markers 2009; 27: 37-44
- Puchau, B; Zulet, MA; Urtiaga, G; Navarro-Blasco, I; Martínez, JA  
**Asymmetric dimethylarginine association with antioxidants intake in healthy young adults: a role as an indicator of metabolic syndrome features**  
Metab. Clin. Exp. 2009; 58: 1483-1488
- Reimann, M; Ziemssen, T; Böger, RH; Schuette, AE  
**Ethnic-specific correlations of visfatin with circulating markers of endothelial inflammation**  
Obesity 2009; 17: 2210-2215.
- Sari, I; Kebapcilar, L; Alacacioglu, A; Bilgir, O; Yildiz, Y; Taylan, A; Yuksel, A; Kozaci, DL  
**Increased levels of asymmetric dimethylarginine (ADMA) in patients with ankylosing spondylitis**  
Intern. Med. 2009; 48: 1363-1368
- Schulze, F; Vollert, J; Maas, R; Müller, R; Schwedhelm, E; Müller, C; Böger, RH; Möckel, M  
**Asymmetric dimethylarginine predicts outcome and time of stay in hospital in patients attending an internal medicine emergency room**  
Clin. Chim. Acta 2009; 401: 20-24
- Sibal, L; Agarwal, SC; Schwedhelm, E; Lüneburg, N; Böger, RH; Home, PD  
**A study of endothelial function and circulating asymmetric dimethylarginine levels in people with Type 1 diabetes without macrovascular disease or microalbuminuria**  
Cardiovasc Diabetol 2009; 8: 27

- Smilde, TD; Damman, K; van der Harst, P; Navis, G; Westenbrink, BD; Voors, AA; Boomsma, F; van Veldhuisen, DJ; Hillege, HL  
**Differential associations between renal function and modifiable risk factors in patients with chronic heart failure**  
Clin Res Cardiol 2009; 98: 121-129
- Tuygun, AK; Keser, M; Tuygun, A; Ketenci, B; Sensöz, Y; Balci, AY; Karaci, AR; Kizilay, M; Yurtseven, N; Yekeler, I  
**Effects of endothelial nitric oxide synthase, interleukin-6 gene polymorphisms and asymmetric dimethylarginine levels on risk factors and lesion sites in peripheral artery disease**  
J. Int. Med. Res. 2009; 37: 1003-1010
- Vladimirova-Kitova, L; Terzieva, D; Marinov, B  
**Intima-media thickness and flow-mediated vasodilation in asymptomatic subjects with newly detected severe hypercholesterolemia**  
Echocardiography 2009; 26: 1060-1068
- Wu, CC; Wen, SC; Yang, CW; Pu, SY; Tsai, KC; Chen, JW  
**Plasma ADMA predicts restenosis of arteriovenous fistula**  
J. Am. Soc. Nephrol. 2009; 20: 213-222
- Young, JM; Terrin, N; Wang, X; Greene, T; Beck, GJ; Kusek, JW; Collins, AJ; Sarnak, MJ; Menon, V  
**Asymmetric dimethylarginine and mortality in stages 3 to 4 chronic kidney disease**  
Clin J Am Soc Nephrol 2009; 4: 1115-1120
- Dimitroulas, T; Giannakoulas, G; Sfetsios, T; Karvounis, H; Dimitroula, H; Koliakos, G; Settas, L  
**Asymmetrical dimethylarginine in systemic sclerosis-related pulmonary arterial hypertension**  
Rheumatology (Oxford). 2008; 47: 1682-1685
- EL-Mesallamy, HO; Abdel Hamid, SG; Gad, MZ  
**Oxidative stress and asymmetric dimethylarginine are associated with cardiovascular complications in hemodialysis patients: improvements by L-arginine intake**  
Kidney Blood Press. Res. 2008; 31: 189-195
- Ferroni, P; Guagnano, MT; Falco, A; Paoletti, V; Manigrasso, MR; Michetti, N; Santilli, F; Guadagni, F; Basili, S; Davì, G  
**Association of low-grade inflammation and platelet activation in patients with hypertension with microalbuminuria**  
Clin. Sci. 2008; 114: 449-455
- Galle, J; Schwedhelm, E; Pinetti, S; Böger, RH; Wanner, CH  
**Antiproteinuric effects of angiotensin receptor blockers: telmisartan versus valsartan in hypertensive patients with type 2 diabetes mellitus and overt nephropathy.**  
Nephrol. Dial. Transplant. 2008; 23: 3174-3183
- Gomes, VA; Casella-Filho, A; Chagas, AC; Tanus-Santos, JE  
**Enhanced concentrations of relevant markers of nitric oxide formation after exercise training in patients with metabolic syndrome**  
Nitric Oxide 2008; 19: 345-350

## 2008

- Alfieri, AB; Briceno, L; Fragasso, G; Spoladore, R; Palloshi, A; Bassanelli, G; Montano, C; Arioli, F; Cuko, A; Ruotolo, G; Margonato, A  
**Differential long-term effects of carvedilol on proinflammatory and antiinflammatory cytokines, asymmetric dimethylarginine, and left ventricular function in patients with heart failure**  
J. Cardiovasc. Pharmacol. 2008; 52: 49-54
- Charitidou, C; Farmakiotis, D; Zournatzi, V; Pidonia, I; Pegiou, T; Karamanis, N; Hatzistilianou, M; Katsikis, I; Panidis, D  
**The administration of estrogens, combined with anti-androgens, has beneficial effects on the hormonal features and asymmetric dimethylarginine levels, in women with the polycystic ovary syndrome.**  
Atherosclerosis 2008; 196: 958-965
- Jacobi, J; Maas, R; Cordasic, N; Koch, K; Schmieder, RE; Böger, RH; Hilgers, KF  
**Role of asymmetric dimethylarginine for angiotensin II-induced target organ damage in mice**  
Am. J. Physiol. Heart Circ. Physiol. 2008; 294: 1058-1066
- Klepárník, M; Pařenica, J  
**HPLC a ELISA stanovení asymetrického dimethylargininu u akutního infarktu myokardu**  
In 52. Studentská vědecká konference - Program a sborník abstrakt. Brno: Lékařská fakulta Masarykovy univerzity 2008; : 34-35
- Korish, AA; Arafah, MM  
**Catechin combined with vitamins C and E ameliorates insulin resistance (IR) and atherosclerotic changes in aged rats with chronic renal failure (CRF).**  
Arch. Gerontol. Geriatr. 2008; 46: 25-39
- Li Volti, G; Sorrenti, V; Acquaviva, R; Murabito, P; Gullo, A; Barcellona, ML; Galvano, F; Rodella, L; Rezzani, R; Vanella, L; Tringali, G; Caruso, M; Gazzolo, D; Di Giacomo, C  
**Effect of ischemia-reperfusion on renal expression and activity of N(G)-N(G)-dimethylarginine dimethylaminohydrolases**  
Anesthesiology 2008; 109: 1054-1062

- Matuszek, MA; Boutcher, SH  
**Elevated levels of circulating cortisol in young normotensive adult men with a family history of hypertension**  
Clin. Exp. Pharmacol. Physiol. 2008; 35: 280-286
  - Onat, A; Hergenç, G; Can, G; Karabulut, A  
**Serum asymmetric dimethylarginine levels among Turks: association with metabolic syndrome in women and tendency to decrease in smokers**  
Turk Kardiyol Dern Ars 2008; 36:
  - Panichi, V; Mantuano, E; Paoletti, S; Santi, S; Manca Rizza, G; Cutrupi, S; Pizzini, P; Spoto, B; Tripepi, G; Zoccali, C  
**Effect of simvastatin on plasma asymmetric dimethylarginine concentration in patients with chronic kidney disease**  
J. Nephrol. 2008; 21: 38-44
  - Potena, L; Fearon, WF; Sydow, K; Holweg, C; Luikart, H; Chin, C; Weisshaar, D; Mocarski, ES; Lewis, DB; Valantine, HA; Cooke, JP  
**Asymmetric dimethylarginine and cardiac allograft vasculopathy progression: modulation by sirolimus**  
Transplantation 2008; 85: 827-833
  - Prefumo, F; Thilaganathan, B; Whitley, GS  
**First-trimester uterine artery resistance and maternal serum concentration of asymmetric dimethylarginine**  
Ultrasound Obstet Gynecol 2008; 31: 153-577
  - Santilli, F; Romano, M; Recchiuti, A; Dragani, A; Falco, A; Lessiani, G; Fioritoni, F; Lattanzio, S; Mattoscio, D; De Cristofaro, R; Rocca, B; Davì, G  
**Circulating endothelial progenitor cells and residual in vivo thromboxane biosynthesis in low-dose aspirin-treated polycythemia vera patients**  
Blood 2008; 112: 1085-1090
  - Sydow, K; Mondon, CE; Schrader, J; Konishi, H; Cooke, JP  
**Dimethylarginine dimethylaminohydrolase overexpression enhances insulin sensitivity**  
Arterioscler. Thromb. Vasc. Biol. 2008; 28: 692-697
  - Teplan, V; Schüeck, O; Racek, J; Lecian, D  
**Asymmetric dimethylarginine in obesity after renal transplantation**  
Journal of Renal Nutrition 2008; 18: 513-520
  - Teplan, V; Schüeck, O; Racek, J; Mareckova, O; Stolova, M; Hanzal, V; Malý, J  
**Reduction of plasma asymmetric dimethylarginine in obese patients with chronic kidney disease after three years of a low-protein diet supplemented with keto-amino acids: a randomized controlled trial**  
Wien. Klin. Wochenschr. 2008; 120: 478-485
  - Teplan, V; Schüeck, O; Racek, J; Siroka, R; Haluzik, M; Kudla, M; Vitko, S  
**Asymmetric dimethylarginine and adiponectin after renal transplantation: role of obesity**  
J Ren Nutr 2008; 18: 154-157
  - Tsukahara, H; Ohta, N; Tokuriki, S; Nishijima, K; Kotsuji, F; Kawakami, H; Ohta, N; Sekine, K; Nagasaka, H; Mayumi, M  
**Determination of asymmetric dimethylarginine, an endogenous nitric oxide synthase inhibitor, in umbilical blood**  
Metab. Clin. Exp. 2008; 57: 215-220
  - Uzun, H; Konukoglu, D; Besler, M; Erdenen, F; Sezgin, C; Muderrisoglu, C  
**The effects of renal replacement therapy on plasma, asymmetric dimethylarginine, nitric oxide and C-reactive protein levels**  
Clin. Invest. Med. 2008; 31: E1-7
  - Vladimirova-Kitova, L; Deneva, T; Angelova, E; Nikolov, F; Marinov, B; Mateva, N  
**Relationship of asymmetric dimethylarginine with flow-mediated dilatation in subjects with newly detected severe hypercholesterolemia**  
Clin Physiol Funct Imaging 2008; 28: 417-425
  - Vostrý, M; Rajdl, D; Eiselt, J; Malánová, L; Píkner, R  
**Adipocytokiny a další prediktory přežití u hemodialyzovaných pacientů s metabolickým syndromem**  
Klin. Biochem. Metab 2008; 16: 239-243
  - Wipff, J; Avouac, J; Borderie, D; Zerkak, D; Lemarechal, H; Kahan, A; Boileau, C; Allanore, Y  
**Disturbed angiogenesis in systemic sclerosis: high levels of soluble endoglin**  
Rheumatology (Oxford) 2008; 47: 972-975
- ## 2007
- 
- Blouet, C; Mariotti, F; Mathe, V; Tome, D; Huneau, JF  
**Nitric oxide bioavailability and not production is first altered during the onset of insulin resistance in sucrose-fed rats**  
Exp. Biol. Med. (Maywood) 2007; 232: 1458-1464
  - CASES, NOND  
**NEW BIOMARKERS IN CORONARY ARTERY DISEASE**  
Cardiovascular Drugs and Therapy 2007; 21: 49-64
  - Devangelio, E; Santilli, F; Formoso, G; Ferroni, P; Bucciarelli, L; Michetti, N; Clissa, C; Ciabattini, G; Consoli, A; Davì, G  
**Soluble RAGE in type 2 diabetes: association with oxidative stress**  
Free Radic. Biol. Med. 2007; 43: 511-518
  - Horowitz, JD; Heresztyn, T  
**An overview of plasma concentrations of asymmetric dimethylarginine (ADMA) in health and disease and in clinical studies: methodological considerations**  
J. Chromatogr. B Analyt. Technol. Biomed. Life Sci. 2007; 851: 42-50
  - Iribarren, C; Husson, G; Sydow, K; Wang, BY; Sidney, S; Cooke, JP  
**Asymmetric dimethyl-arginine and coronary artery calcification in young adults entering middle age: the CARDIA Study**  
Eur J Cardiovasc Prev Rehabil 2007; 14: 222-229

- Konishi, H; Sydow, K; Cooke, JP  
**Dimethylarginine dimethylaminohydrolase promotes endothelial repair after vascular injury**  
J. Am. Coll. Cardiol. 2007; 49: 1099-1105
- Konukoglu, D; Uzun, H; Firtina, S; Cigdem Arica, P; Kocael, A; Taskin, M  
**Plasma adhesion and inflammation markers: asymmetrical dimethyl-L-arginine and secretory phospholipase A2 concentrations before and after laparoscopic gastric banding in morbidly obese patients**  
Obes Surg. 2007; 17: 672-678
- Melikian, N; Wheatcroft, SB; Ogah, OS; Murphy, C; Chowiecnyk, PJ; Wierzbicki, AS; Sanders, TA; Jiang, B; Duncan, ER; Shah, AM; Kearney, MT  
**Asymmetric dimethylarginine and reduced nitric oxide bioavailability in young Black African men**  
Hypertension 2007; 49: 873-877
- Öguz, A; Uzunlulu, M; Yorulmaz, E; Yalcin, Y; Hekim, N; Fici, F  
**Effect of nebivolol and metoprolol treatments on serum asymmetric dimethylarginine levels in hypertensive patients with type 2 diabetes mellitus**  
Anadolu Kardiyol Derg 2007; 7: 383-7
- Santilli, F; Bucciarelli, L; Noto, D; Cefalù, AB; Davì, V; Ferrante, E; Pettinella, C; Aversa, MR; Ciabattone, G; Davì, G  
**Decreased plasma soluble RAGE in patients with hypercholesterolemia: effects of statins**  
Free Radic. Biol. Med. 2007; 43: 1255-1262
- Shankar, SS; Considine, RV; Clauss, M  
**Effect of Free Fatty Acids on Plasma and Endothelial Cell ADMA Levels**  
Diabetes Jun 2007; 56: 191
- Siroka, R; Trefil, L; Rajdl, D; Racek, J; Cibulka, R  
**Asymmetric dimethylarginine – comparison HPLC and ELISA method (Letter to the Editor)**  
J. Chromatogr. 2007; B850: 586-587
- Soveri, I; Lind, L; Wikstrom, B; Zilmer, M; Zilmer, K; Fellström, B  
**Improvement in central arterial pressure waveform during hemodialysis is related to a reduction in asymmetric dimethylarginine (ADMA) levels**  
Nephron Clin Pract 2007; 106: 180-186
- Spoto, B; Parlongo, RM; Parlongo, G; Sgro', E; Zoccali, C  
**The enzymatic machinery for ADMA synthesis and degradation is fully expressed in human adipocytes**  
J. Nephrol. 2007; 20: 554-559
- Thum, T; Fleissner, F; Klink, I; Tsikas, D; Jakob, M; Bauersachs, J; Stichtenoth, DO  
**Growth hormone treatment improves markers of systemic nitric oxide bioavailability via insulin-like growth factor-I**  
J. Clin. Endocrinol. Metab. 2007; 92: 4172-4179
- Tousoulis, D; Antoniades, C; Vasiliadou, C; Kourtellis, P; Koniari, K; Marinou, K; Charakida, M; Ntarladimas, I; Siasos, G; Stefanadis, C  
**Effects of atorvastatin and vitamin C on forearm hyperaemic blood flow, asymmetrical dimethylarginine levels and the inflammatory process in patients with type 2 diabetes mellitus**  
Heart 2007; 93: 244-246
- Tsukahara, H; Ohta, N; Tokuriki, S; Nishijima, K; Kotsuji, F; Kawakami, H; Ohta, N; Sekine, K; Nagasaka, H; Mayumi, M  
**Determination of asymmetric dimethylarginine, an endogenous nitric oxide synthase inhibitor, in umbilical blood**  
Metabolism Clin. Exp. 2007; 57: 215-220
- Vizzutti, F; Romanelli, RG; Arena, U; Rega, L; Brogi, M; Calabresi, C; Masini, E; Tarquini, R; Zipoli, M; Boddi, V; Marra, F; Laffi, G; Pinzani, M  
**ADMA correlates with portal pressure in patients with compensated cirrhosis**  
Eur. J. Clin. Invest. 2007; 37: 509-15
- Vladimirova-Kitova, L; Terzieva, D; Nikolov, F  
**Asymmetric dimethylarginine in hypercholesterolemic patients**  
Journal of IMAB 2007; 13: 110-112
- Vladimirova-Kitova, L; Terzieva, D; Nikolov, F; Mateva, N  
**APOLIPOPROTEIN-B AS A PREDICTOR OF ASYMMETRIC DIMETHYLARGININE IN HYPERCHOLESTEROLEMIC PATIENTS**  
J of IMAB 2007; 13: 113-115
- Wanby, P; Nilsson, I; Brudin, L; Nyhammar, I; Gustafsson, I; Carlsson, M  
**Increased plasma levels of asymmetric dimethylarginine in patients with carotid stenosis: no evidence for the role of the common FABP2 A54T gene polymorphism**  
Acta Neurol. Scand. 2007; 115: 90-96
- Weber, T; Maas, R; Auer, J; Lamm, G; Lassnig, E; Rammer, M; O'Rourke, MF; Böger, RH; Eber, B  
**Arterial wave reflections and determinants of endothelial function a hypothesis based on peripheral mode of action**  
Am. J. Hypertens. 2007; 20: 256-262
- Xu, L; Wang, B; Kaur, K; Kho, MF; Cooke, JP; Giffard, RG  
**NOx and ADMA changes with focal ischemia, amelioration with the chaperonin GroEL**  
Neurosci. Lett. 2007; 418: 201-4

## 2006

- Antoniades, C; Tousoulis, D; Marinou, K; Vasiliadou, C; Tentolouris, C; Bouras, G; Pitsavos, C; Stefanadis, C  
**Asymmetrical dimethylarginine regulates endothelial function in methionine-induced but not in chronic homocystinemia in humans: effect of oxidative stress and proinflammatory cytokines**  
Am. J. Clin. Nutr. 2006; 84: 781-788

- Dooley, A; Gao, B; Bradley, N; Abraham, DJ; Black, CM; Jacobs, M; Bruckdorfer, KR  
**Abnormal nitric oxide metabolism in systemic sclerosis: increased levels of nitrated proteins and asymmetric dimethylarginine**  
Rheumatology (Oxford) 2006; 45: 676-684
  - El-Saleh, SC  
**Protection by Natural Honey Against Hyperhomocysteinemia in Rats**  
Vascular Disease Prevention 2006; 3: 313-318
  - Jacobi, J; Porst, M; Cordasic, N; Namer, B; Schmieder, RE; Eckardt, KU; Hilgers, KF  
**Subtotal nephrectomy impairs ischemia-induced angiogenesis and hindlimb re-perfusion in rats**  
Kidney Int. 2006; 69: 2013-2021
  - O'Dwyer, MJ; Dempsey, F; Crowley, V; Kelleher, DP; McManus, R; Ryan, T  
**Septic shock is correlated with asymmetrical dimethyl arginine levels, which may be influenced by a polymorphism in the dimethylarginine dimethylaminohydrolase II gene: a prospective observational study**  
Crit Care 2006; 10: 139
  - Reimann, M; Dierkes, J; Carlsohn, A; Talbot, D; Ferrari, M; Hallund, J; Hall, WL; Vafeiadou, K; Huebner, U; Branca, F; Bugel, S; Williams, CM; Zunft, HJ; Koebnick, C  
**Consumption of soy isoflavones does not affect plasma total homocysteine or asymmetric dimethylarginine concentrations in healthy postmenopausal women**  
J. Nutr. 2006; 136: 100-105
  - Reimann, M; Vafeiadou, K; Hall, WL; Dierkes, J; Nilsson, M; Dahlman-Wright, K; Ferrari, M; Huebner, U; Hallund, J; Bugel, S; Branca, F; Williams, CM; Koebnick, C  
**Evidence for associations between common polymorphisms of estrogen receptor beta gene with homocysteine and nitric oxide**  
Climacteric 2006; 9: 215-223
  - Schulze, F; Lenzen, H; Hanefeld, C; Bartling, A; Osterziel, KJ; Goudeva, L; Schmidt-Lucke, C; Kusus, M; Maas, R; Schwedhelm, E; Strödter, D; Simon, BC; Mügge, A; Daniel, WG; Tillmanns, H; Maisch, B; Streichert, T; Böger, RH  
**Asymmetric dimethylarginine is an independent risk factor for coronary heart disease: results from the multicenter Coronary Artery Risk Determination Investigating the Influence of ADMA Concentration (CARDIAC) study**  
Am. Heart J. 2006; 152: 493-498
  - Soveri, I; Holdaas, H; Jardine, A; Gimpelewicz, C; Staffler, B; Fellström, B  
**Renal transplant dysfunction--importance quantified in comparison with traditional risk factors for cardiovascular disease and mortality**  
Nephrol. Dial. Transplant. 2006; 21: 2282-2290
  - Wang, TD; Chen, WJ; Cheng, WC; Lin, JW; Chen, MF; Lee, YT  
**Relation of improvement in endothelium-dependent flow-mediated vasodilation after rosiglitazone to changes in asymmetric dimethylarginine, endothelin-1, and C-reactive protein in nondiabetic patients with the metabolic syndrome**  
Am. J. Cardiol. 2006; 98: 1057-1062
  - Wierzbicki, AS; Solomon, H; Lumb, PJ; Lyttle, K; Lambert-Hammill, M; Jackson, G  
**Asymmetric dimethyl arginine levels correlate with cardiovascular risk factors in patients with erectile dysfunction**  
Atherosclerosis 2006; 185: 421-425
- ## 2005
- 
- Jacobi, J; Sydow, K; von Degenfeld, G; Zhang, Y; Dayoub, H; Wang, B; Patterson, AJ; Kimoto, M; Blau, HM; Cooke, JP  
**Overexpression of dimethylarginine dimethylaminohydrolase reduces tissue asymmetric dimethylarginine levels and enhances angiogenesis**  
Circulation 2005; 111: 1431-1438
  - Krempf, TK; Maas, R; Sydow, K; Meinertz, T; Böger, RH; Kähler, J  
**Elevation of asymmetric dimethylarginine in patients with unstable angina and recurrent cardiovascular events**  
Eur. Heart J. 2005; 26: 1846-1851
  - Martens-Lobenhoffer, J; Westphal, S; Awiszus, F; Bode-Böger, SM; Luley, C  
**Determination of asymmetric dimethylarginine: liquid chromatography-mass spectrometry or ELISA?**  
Clin. Chem. 2005; 51: 2188-2190
  - Ravani, P; Tripepi, G; Malberti, F; Testa, S; Mallamaci, F; Zoccali, C  
**Asymmetrical dimethylarginine predicts progression to dialysis and death in patients with chronic kidney disease: a competing risks modeling approach**  
J. Am. Soc. Nephrol. 2005; 16: 2449-2455
  - Schnabel, R; Blankenberg, S; Lubos, E; Lackner, KJ; Rupprecht, HJ; Espinola-Klein, C; Jachmann, N; Post, F; Peetz, D; Bickel, C; Cambien, F; Tiret, L; Münzel, T  
**Asymmetric dimethylarginine and the risk of cardiovascular events and death in patients with coronary artery disease: results from the AtheroGene Study**  
Circ. Res. 2005; 97: 53-59
  - Schulze, F; Maas, R; Freese, R; Schwedhelm, E; Silberhorn, L; Böger, RH  
**Determination of a reference value for N,N-dimethyl-L-arginine in 500 subjects**  
Eur. J. Clin. Invest. 2005; 35 : 622-626



- Šíroká, R; Racek, J; Filipovský, J  
**Asymetrický dimethylarginin (ADMA)**  
Klin Biochem Metab 2005; 13: 131-134
- Šíroká, R; Trefil, L; Rajdl, D; Racek, J; Rusňáková, H  
**Asymetrický dimethylarginin, homocystein a funkce ledvin**  
Klin. Biochem. Metab. 2005; 13: 184-186
- Tanaka, M; Sydow, K; Gunawan, F; Jacobi, J; Tsao, PS; Robbins, RC; Cooke, JP  
**Dimethylarginine dimethylaminohydrolase overexpression suppresses graft coronary artery disease**  
Circulation 2005; 112: 1549-1556
- Valtonen, P; Karppi, J; Nyssönen, K; Valkonen, VP; Halonen, T; Punnonen, K  
**Comparison of HPLC method and commercial ELISA assay for asymmetric dimethylarginine (ADMA) determination in human serum**  
J. Chromatogr. B Analyt. Technol. Biomed. Life Sci. 2005; 828: 97-102

## 2004

---

- Schulze, F; Wesemann, R; Schwedhelm, E; Sydow, K; Albsmeier, J; Cooke, JP; Böger, RH  
**Determination of ADMA using a novel ELISA assay**  
Clin. Chem. Lab. Med. 2004; 42: 1377-1383
- Trapp, G; Sydow, K; Dulay, MT; Chou, T; Cooke, JP; Zare, RN  
**Capillary electrophoretic and micellar electrokinetic separations of asymmetric dimethyl-L-arginine and structurally related amino acids: quantitation in human plasma**  
J Sep Sci 2004; 27: 1483-1490