

1 COMPARISON STUDIES

Mediagnost Adiponectin, E09 was compared with two different, commercially available test systems. In linear regression analysis both tests showed a good coefficient of determination ($R^2=0.896$ and $R^2=0.97$). Thus, the comparability of the results between the Mediagnost E09 and other test systems is possible. In dependence on the respective system the absolute deviation of measurement results is different, but because of the excellent correlation the results are comparable after applying a factor. The results of both studies are shown in figure 1.

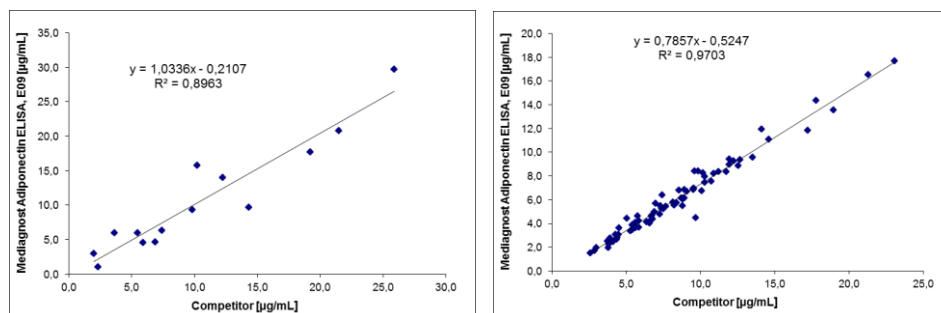


Figure 1 Comparison of Mediagnost E09 with commercially available test systems (left) radioimmunoassay (n=14) LINCO and (right) **ELISA R&D** (n=84)

Samples suitable for scientific application

Serum, plasma, saliva, urine, breast milk and cell culture supernatant of human cell Lines. The recommended dilution for serum and plasma samples in Dilution Buffer **VP: (1:310)**. In the other samples, the Adiponectin levels can vary considerable, the optimal dilution must be found out by the customer.

Table 7 Results of sample matrix tests . Adiponectin was added to the respectively diluted samples. Enriched samples were measured without further dilution. Shown is the relative recovery of added Adiponectin of the concentration measured in respectively enriched assay buffer.

Matrix Dilution	Urine	Saliva	Breast Milk	Cell Culture Medium 10% FCS	Cell Culture Medium
1:2	80	87	89	83	95
1:5	95	80	92	92	97
1:10	92	87	-	101	85
1:20	94	99	-	83	91