



Antioxidant Capacity Potential Assay

Catalog Number: PAC39-K01
96 Wells
For Research Use Only
v. 1.0

EAGLE BIOSCIENCES, INC.
20A NW Blvd, Suite 112
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



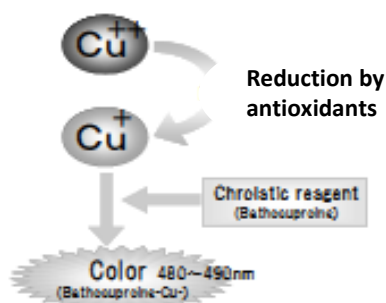
Intended Use:

The Eagle Biosciences Antioxidant Capacity Potential Assay kit is intended for the quantitative determination of antioxidant capacity in biological samples as well as well food and beverage samples by enzyme linked immunoassay (ELISA). The Antioxidant Capacity Potential Assay kit is for research use only and not to be used in diagnostic procedures.

Assay Background:

Oxidative stress plays an important role in various diseases and aging. The control of oxidative stress is expected to be useful to prevent diseases and aging. Oxidative stress is caused by the imbalance between reactive oxygen species (ROS) and antioxidant defense system. For accurate assessment of oxidative stress, measurement of ROS, oxidative damage and antioxidant activity may be essential. The Eagle Biosciences Potential Antioxidant Capacity Assay can detect not only hydrophilic antioxidants such as Vitamin C and glutathione, but also can detect hydrophobic antioxidants such as Vitamin E. This Antioxidant Capacity Potential Assay kit is applicable for assessment of total antioxidants of serum, foods and beverage samples.

Principle of Procedure:



Samples are mixed with Cu^{++} Solution. Cu^{++} are reduced by antioxidants to form Cu^{+} . Reduced Cu^{+} react with Chromatic Solution (Bathocuproine), and can be detected by absorbance at wavelength 480 to 490 nm. Antioxidant capacity can be calculated from the Cu^{+} formed.

Materials Provided with the Assay:

- | | | |
|-------------------------------|------------------|--------------------------------|
| • Standard (Uric acid powder) | 1 vial (for 2mM) | Dissolve with distilled water. |
| • Sample diluent | 60 mL x 1 bottle | Ready to use. |
| • Cu^{++} solution | 5 mL x 1 bottle | Ready to use. |
| • Stop solution | 5 mL x 1 bottle | Ready to use. |
| • Micro titer plate | 1 plate | |



Materials Required but not Provided:

- Microplate reader (measuring wavelength 490 nm)
- Pipettes and pipette tips
- Plastic test tubes
- Distilled water
- NaOH, HCl solution and pH meter (Not required if standards are prepared with distilled water only)

Specifications:

- Assay range: 21.9~4378 $\mu\text{mol/L}$ (Cupric ion reducing power)
- Storage of the Antioxidant Capacity Potential Assay: Room Temperature
- Expiration date: 3 years from manufacture date

Assay Procedure

1. Reconstitute of Standard (2mM Uric acid solution).
 - There are two options for preparation. Select one.
 - i. Option 1: Add distilled water to the Standard vial, and stand for 3 or 4 hours at room temperature. The volume of distilled water is indicated on the label of the vial.
 - ii. Option 2: If you wish to prepare standard solution immediately, add 1mL of 10% (w/v) NaOH to Standard vial and dissolve completely, followed by pH adjustment (pH7.4) by HCl solution. Add distilled water to make the total volume as indicated on the label. 2mM uric acid solution can be stored at below -70°C for 1 year.
2. Preparation of standards.
 - Dilute 2mM uric acid solution with distilled water for 2, 4, 8, 16 and 32 times, result in 5 levels of diluted standards (1 mM, 0.5 mM, 0.25 mM, 0.125 mM and 0.063 mM respectively).
3. Preparation of samples.
 - For serum samples, fresh frozen samples are recommended because some antioxidants such as vitamin C, uric acid and coenzyme Q10 are unstable. For other samples types such as beverages and food, see section "Assay Examples", and dilute with distilled water.



4. Assay procedure.

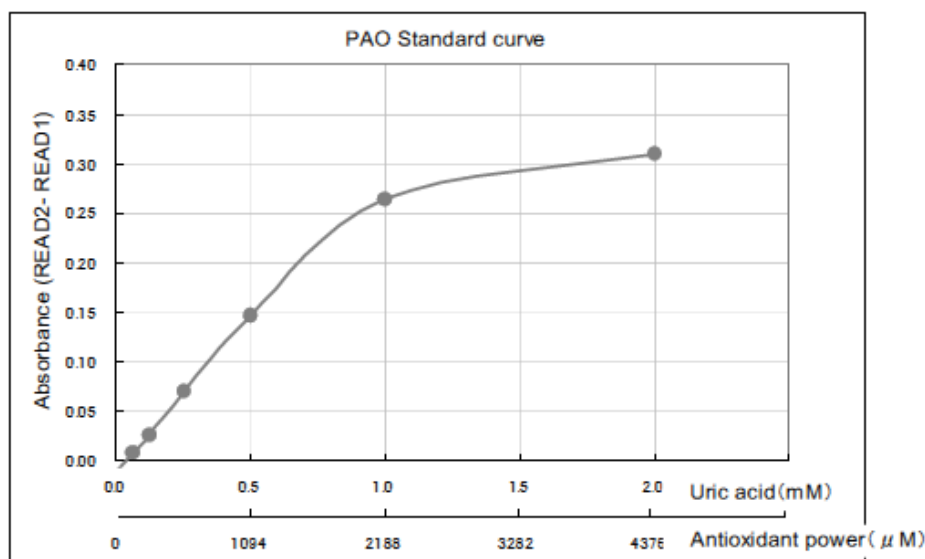
1. Please prepare plastic test tubes for 6 levels of standards and each sample. Pour 390 μL of Sample Diluent, and add 10 μL of standards or diluted samples.
2. Pour 200 μL of mixture to Micro titer plate. Use 200 μL of Sample Diluent for blank well.
3. Read absorbance at 490 nm (as READ1).
4. Add 50 μL of Cu^{++} solution to each well, mix gently, and incubate at room temperature for 3 minutes.
5. Add 50 μL of Stop solution, mix gently, and read absorbance at 490 nm (as READ2).

5. Determination of antioxidant power of samples.

- Please draw standard curves by plotting the difference of absorbance readings ($\text{READ2} - \text{READ1}$) as vertical axis, and concentration of uric acid standards (mM) as horizontal axis.
- Calculate the corresponding uric acid concentration of samples. Multiply corresponding uric acid concentration (mM) of samples by 2189, to estimate antioxidant power ($\mu\text{mol/L}$).

1mM of uric acid = 2189 $\mu\text{mol/L}$ (copper reducing power)

Typical Standard Curve





Assay Examples:

Sample	Pre-dilution	Antioxidant power (μmol/L)
Human serum	Not required	1069±145
Human urine	Mix with 3 volumes of Distilled Water	5508
Red wine	Mix with 7 volumes of Distilled Water	45479
Japanese SAKE (rice wine)	Not required	18~211
Black tea	Mix with 7 volumes of Distilled Water	TBD
Coffee	Mix with 27 volumes of Distilled Water	TBD
Green tea	Mix with 7 volumes of Distilled Water	8728~46687

Additional dilution is recommended if the antioxidant power is over 2000 μmol/L antioxidant power. For example, some green tea products which contain high concentration of catechin should be diluted by 40 times (mix 1 volume of sample and 39 volume of distilled water). Some samples which contain chelating agents such as EDTA can't be applied with this Antioxidant Capacity Potential Assay.

References:

- Oxidative stress and its association with coronary artery disease and different atherogenic risk factors C. VASSALLE, L. PETROZZI, N. BOTTO, M. G. ANDREASSI & G. C. ZUCHELLI Journal of Internal Medicine 256, p308–315 (2004)
- Oxidative imbalance and cathepsin D changes as peripheral blood biomarkers of Alzheimer disease: A pilot study E Strafacea, P Matarrese, L Gambardella, R Vona, A Sgadari, MC Silveri, W Malorni FEBS Letters 579, p2759-766 (2005)
- Restored Antioxidant Capacity Parallels the Immunologic and Virologic Improvement in Children with Perinatal Human Immunodeficiency Virus Infection Receiving Highly Active Antiretroviral Therapy.
- M Martino, F Chiarelli, M Moriondo, M Torello, C Azzari, and L Galli. Clinical Immunology, Vol.100(1),p82-6 (2001)
- Antioxidant capacity as a reliable marker of stress in dairy calves transported by road. P Preget, E Bollo, FT Cannizzo, B Biolatti, E Contato, PG Biolatti: Veterinary Record Vol. 156, p53-54 (2005)

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This



warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.