

OxiSelect™ Total Antioxidant Capacity (TAC) Assay Kit

The OxiSelect™ Total Antioxidant Capacity (TAC) Assay measures the total antioxidant capacity of biomolecules from a variety of samples via a SET mechanism. In the presence of antioxidants, copper(II) is reduced to copper(I). In turn, the copper(I) ions react with a chromogen to produce a color with maximum absorbance at 490nm.

Ordering Information

[Order Online »](#)

CBI-STA-360	200 assays
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Measures total antioxidant capacity based on reduction of copper(II) to copper(I)
- Suitable for use with plasma, serum, urine, cell lysates, tissue homogenates and food extracts
- Works with a wide variety of antioxidants

Handling & Storage

Use/Stability Upon receipt store the Reaction Buffer (100X) at 4°C. Store all remaining kit components at room temperature.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)