

OxiSelect™ Ascorbic Acid Assay Kit (FRASC)

Ascorbic acid is a vital water-soluble antioxidant found in living organisms. Ascorbic acid is critical for a variety of functions related to tissue growth and wound healing, neurotransmitter formation, blood cholesterol levels, as well as free radical neutralization. The OxiSelect™ Ascorbic Acid Assay Kit is a quantitative assay for measuring ascorbic acid within various samples. The assay is based on the Ferric Reducing/Antioxidant Ascorbic Acid (FRASC) chemistry driven by the electron donating reducing power of antioxidants. The kit employs ascorbate oxidase, which allows the user to differentiate the ascorbic acid content from other antioxidants present within their sample. Ascorbic acid levels in a sample are determined by measuring the difference in optical density between two sample wells, one with and one without the enzyme.

Ordering Information

[Order Online »](#)

CBI-STA-860	200 assays
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Specifically differentiates ascorbic acid content from other antioxidant compounds
- Detects ascorbic acid concentrations as low as 1 μ M
- Suitable for use with serum, plasma, urine, saliva, tissue homogenates, cell extracts, and purified food or drug extracts
- NOTE: Each sample replicate requires 2 assays for proper data calculation.

Handling & Storage

Use/Stability Upon receipt, prepare single use aliquots and store the Ascorbate Oxidase at -80°C. Store the remaining kit components at 4°C.

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)