

Protein Carbonyl Assay

A plate based colorimetric assay to measure protein carbonyls in a variety of samples with a sensitivity well below 1 nmole.

Oxidative stress due to an imbalance between ROS production and antioxidant defense leads to cellular damage including the oxidation of amino acid side chains, resulting in protein carbonyls. Protein carbonylation disrupts the function and structure of proteins involved in various metabolic pathways. Protein carbonylation triggers immune responses, leading to chronic inflammation affecting things like insulin sensitivity and lipid metabolism. Elevated levels of protein carbonyls are associated with age-related diseases like neurodegenerative disorders, cardiovascular disease, and diabetes. AkrivisBio's Protein Carbonyl Assay is a simple sensitive means of determining protein carbonyls in a variety of samples with a sensitivity well below 1 nmole.

Ordering Information

[Order Online »](#)

AKR-MA-0150	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Species Reactivity Species independent

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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