

# Protein carbonyl ELISA kit

**High sensitivity ELISA for quantitative determination of carbonylated protein levels in plasma, other body fluids, cell and tissue extracts.**

The Protein carbonyl ELISA kit is an antibody-based detection set. It has higher sensitivity, lower background, is less labour-intensive and handles more samples per day as compared to colorimetric assays.

Protein carbonyls are formed by a variety of oxidative mechanisms and are sensitive indices of oxidative injury. The quantity of protein carbonyls in a protein sample can be determined by derivatising with dinitrophenylhydrazine (DNP) and measure bound DNP colorimetrically or immunologically. The ELISA method enables carbonyls to be measured quantitatively with microgram quantities of protein.

Citations: 44

[View Online »](#)

## Ordering Information

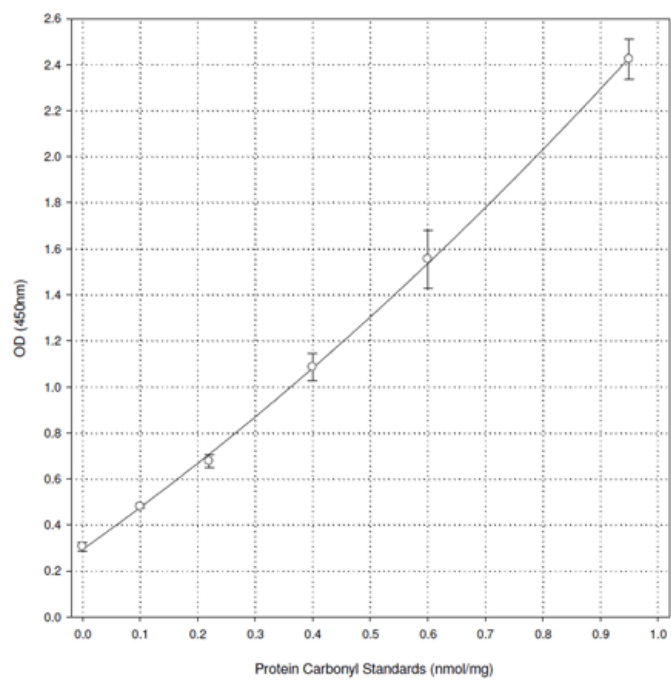
[Order Online »](#)

ALX-850-312-KI01	96 wells
------------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- High sensitivity ELISA, detecting low levels of carbonylated proteins
- Validated in various sample types
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Standard curve of Protein carbonyl ELISA kit

## Handling & Storage

Use/Stability	Remove box containing components G, K, and L and store at -20°C! Remove box containing components H and I and store at 4°C! DO NOT FREEZE! Store the remaining of the kit at room temperature. The kit is stable at room temperature for up to 6 weeks.
Handling	Do not freeze.
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	ELISA
Application Notes	For quantitative determination of carbonylated protein levels in plasma, other body fluids, cell and tissue extracts.
Assay Time	4.5 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Quantity	96 wells (~80 tests)
Species Reactivity	Species independent
Wavelength	450 nm

Last modified: May 29, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)