

Cu/Zn-Superoxide Dismutase ELISA kit

Late stages of apoptosis are characterized by the loss of membrane integrity and a release of the intracellular Cu/Zn SOD.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-0330	96 wells
--------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	SOD (Cu/Zn)
Application	ELISA
Application Notes	For the quantitative determination of human Cu/Zn SOD in cell culture supernatants, serum and other body fluids.
Assay Time	<2 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Quantity	96 wells (~80 tests).
Sensitivity	0.04ng/ml (range 0.08 to 5ng/ml)
Species Reactivity	Human
UniProt ID	P00441
Wavelength	450 nm

Last modified: May 29, 2024



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