

Ceruloplasmin (human plasma)

Serum copper transport protein

Serum copper transport and iron-oxidizing protein. Expressed in plasma at concentrations of 1-5 μ M. Contains six copper centers and is known to oxidize amines in a process coupled to the reduction of molecular oxygen. Has ferroxidase activity that is responsible for the oxidation of ferrous iron to its ferric form, which is necessary for efficient iron efflux from the cell (e.g. during hypoxia). Has previously been considered a target for nitric oxide (NO) and an inhibitor of endothelial nitric oxide synthase (eNOS/NOS III). Catalyzes S-nitrosothiol formation in cell culture media. Is a NO oxidase and nitrite synthase that determines endocrine NO homeostasis.

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-089-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ferroxidase

CAS 9031-37-2

Formulation Lyophilized from 50mM potassium phosphate, pH 6.8, containing 100mM potassium chloride, 20mM ϵ -aminocaproic acid and 5mM EDTA.

MI 14: 2006

MW ~134kDa

Purity $\geq 95\%$ (SDS-PAGE)

Purity Detail Plasma has been tested negative for HBsAg and for antibodies to HIV and HCV.

Reconstitution Reconstitute with distilled water.

Source Isolated from human plasma.

UniProt ID P00450



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)