

Myeloperoxidase (human leukocytes)

Haem protein that catalyzes oxidations by H_2O_2 , including MPO-chloride-mediated killing of microbes and tumor cells, inactivation of chemotactic factors, cross-linking of proteins and iodination of proteins.

Ordering Information

[Order Online »](#)

ALX-200-600A-U025	25µg
ALX-200-600A-5025	5x25µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stored at -20°C this product is stable for years. Storage of the reconstituted form at 4°C for 2 weeks has no effect on enzyme activity. If solutions must be stored for extended periods of time, protein concentration should be kept above 1mg/ml and the activity should be determined prior to use. Freezing solutions of Myeloperoxidase may cause substantial loss of activity.
Short Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	MPO
Appearance	Green solid
CAS	9003-99-0
Formulation	Lyophilized from 5mM phosphate buffer, pH 7.0.
Purity	≥98%
Reconstitution	Reconstitution of 25 µg with 100 µl distilled water gives approximately 3.45µM calculated with a molar extinction coefficient of 91.000 M-1cm-1. per hem at 429nm.
Source	Isolated from human leukocytes. Negative for HBsAG-, HCV- and HIV-antibodies.
UniProt ID	P05164



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)