

Blocking peptide for Heme oxygenase 1 pAb (Prod. No. BML- HC3001)

Synthetic peptide corresponding to amino acid residues [12-25] + Cys-NH₂ of human heme oxygenase-1. Complementary peptide to rabbit polyclonal antibody (BML-HC3001) for use in control experiments. Preincubation of antibody with this peptide abolishes the antibody reactivity with HO-1 protein.

Ordering Information

[Order Online »](#)

BML-HP9301-0100	100µg
-----------------	-------

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

Appearance Lyophilized solid.

MW 1671.9

Purity ≥96%

Sequence H-Asp-Leu-Ser-Glu-Ala-Leu-Lys-Glu-Ala-Thr-Lys-Glu-Val-His-Cys-NH₂

Solubility Soluble in Water.

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)