

Hemopexin (human) monoclonal antibody (4) (biotin conjugate)

Hemopexin is a serum glycoprotein that binds heme and transports it to the liver for breakdown and iron recovery, after which the free hemopexin returns to the circulation. Structurally, hemopexin consists of two similar halves of approximately 200 amino acid residues connected by a histidine-rich hinge region. Each half is formed by the repetition of a basic unit of some 35 to 45 residues.

This antibody is covered by our [Worry-Free Guarantee](#).

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	HPX, Pexin, β -1B-glycoprotein
Application	ELISA
Clone	4
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
GenBank ID	3263
Host	Mouse
Immunogen	Purified human hemopexin from plasma.
Isotype	IgG2bk
Purity Detail	Protein A/G-purified.
Species Reactivity	Human
UniProt ID	P02790
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

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

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)