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 histidinerich 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.

Ordering Information

Order Online »

BPD-ABS-013-04B-015

150µg

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

