HO-1 monoclonal antibody (HO-1-2)

Heme Oxygenase-1 (HO-1) also known as Hsp32, is the inducible isoform of heme oxygenase that catalyzes the NADPH, oxygen, and cytochrome P450 reductase dependent oxidation of heme to carbon monoxide, ferrous iron and biliverdin which is rapidly reduced to bilirubin. These products of the HO reaction have important physiological effects: carbon monoxide is a potent vasodilator and has been implicated to be a physiological regulator of cGMP and vascular tone; biliverdin and its product bilirubin are potent antioxidants; "free" iron increases oxidative stress and regulates the expression of many mRNAs (e.g., DCT-1, ferritin and transferring receptor) by affecting the conformation of iron regulatory protein (IRP)-1 and its binding to iron regulatory elements (IREs) in the 5'- or 3'- UTRs of the mRNAs. To date, three identified heme oxygenase isoforms are part of the HO system that catalyze heme into biliverdin and carbon monoxide. These are inducible HO-1 or Hsp32, constitutive HO-2 that is abundant in the brain and testis, and HO-3 which is related to HO-2 but is the product of a different gene. The HO system is the rate-limiting step in heme degradation and HO activity decreases the levels of heme which is a well known potent catalyst of lipid peroxidation and oxygen radical formation.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

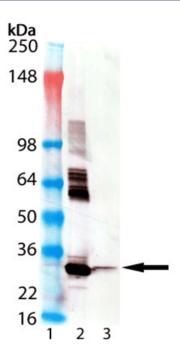
Order Online »

ENZ-ABS687-0200

200µg

Manuals, SDS & CofA

View Online »



Western blot analysis of ENZ-ABS687: Lane 1: MW Marker, Lane 2: HO-1 Recombinant Rat Protein (Prod. No. ADI-SPP-730), Lane 3: Rat Liver Microsomes (Prod. No. ADI-LYT-RM100).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HMOX1, Hsp32, Heat shock protein 32, Heme oxygenase

1

Application Flow Cytometry, IHC, WB

Application NotesDetects a band of ~32kDa by Western blot.

Clone HO-1-2

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

GenBank ID J02722

Host Mouse

Immunogen Native rat HO-1.

lgG2b

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:100)Western Blot (1:1,000)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Dog, Guinea pig, Hamster, Human, Monkey, Mouse,

Porcine, Rabbit, Rat

UniProt ID P06762

Last modified: May 29, 2024



info-

eu@enzolifesciences.com