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

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.

Citations: 2

View Online »

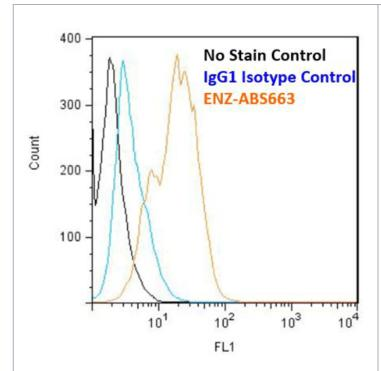
Ordering Information

Order Online »

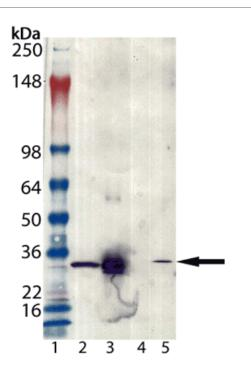
ENZ-ABS663-1000	1mg
ENZ-ABS663-0050	50µg
ENZ-ABS663-0200	200μg

Manuals, SDS & CofA

View Online »



Flow cytometry analysis of 10^6 Jurkat cells stained using HO-1 monoclonal antibody (HO-1-1), (Prod. No. ENZ-ABS663) conjugated to FITC, at a concentration of $10\mu g/ml$.



Western Blot analysis of ENZ-ABS663: Lane 1: MW Marker, Lane 2: HO-1 (rat), (recombinant) (Prod. No. ADI-SPP-730), Lane 3: HO-1 (human), (recombinant) (Prod. No. ADI-SPP-732), Lane 4: HO-2 (human), (recombinant) (Prod. No. ADI-SPP-550), Lane 5: MDBK cell lysate.

Handling & Storage

Handling Avoid freeze/thaw cycles.

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, ICC, IHC (PS), WB

Clone HO-1-1

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

sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to the sequence near the

N-terminus of human HO-1.

Isotype IgG1ĸ

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (10µg/ml)Western Blot (1:1,000

ECL)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 Bovine, Dog, Human, Mouse, Rat

UniProt ID P09601

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com