Heme oxygenase 2 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

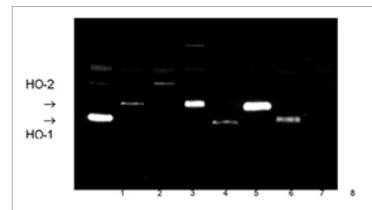
Ordering Information

Order Online »

BML-HC3002-0025	25µl
BML-HC3002-0100	100µl

Manuals, SDS & CofA

View Online »



Western blot analysis: HO-1 was detected using HO-1, pAb (Prod. no. BML-HC3001) (lanes 1, 3, 5 and 7), and HO-2 was detected using HO-2, pAb (Prod. no. BML-HC3002) (lanes 2, 4, 6 and 8). Both antibodies were used at a dilution of 1:1000. Luminograph (ECL, 15 sec. exposure, ref. FM020/054) showing HO-1 and HO-2 expression in lysates prepared from gIFN/LPS stimulated RAW264.7 mouse macrophage cells (lanes 1 and 2), HepG2 (lanes 3 and 4), rat testis (HP9312, lanes 5 and 6) and rat spleen (HP9313, lanes 7 and 8).

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 HMOX2, Hsp32, Heat shock protein 32, HO-2

Application IHC, WB

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 246-264 of human

heme oxygenase 2 (HO-2) protein (E.C. 1.14.99.3).

Purity Detail Protein G-affinity purified.

Source Purified from rabbit serum.

Species Reactivity Human, Rat

Technical Info / Product Notes

The antiserum was raised to a synthetic peptide, sequence: H-E-T-L-E-D-G-F-P-V-H-D-G-K-G-D-M-R-K-C-NH₂ corresponding to amino acid residues 246-264, in amidated form, of the human HO-2 protein (McCoubrey et al., 1992). Two amino acid substitutions (Phe252 -> Leu; Met261 -> Val) exist in the aligned region within rat HO-2, and four (Leu248 -> Val; Gly251 -> Arg; Phe252 -> Leu; Met261 -> Val) in rabbit HO-2. The peptide was conjugated at the C-terminal cysteine residue via a maleimido linkage to keyhole limpet haemocyanin. The antiserum has been purified by caprylic acid and ammonium sulphate precipitation, and by protein G affinity chromatography.

Application: Heme oxygenase-2 is constitutively expressed in many tissues, but the protein levels are extremely variable. The use of enriched enzyme preparations is recommended for experimental analysis. This antibody has been shown to react strongly with HO-2 on Western blots of microsomal preparations of unstimulated rat brain (HP 9311), testis (HP 9312) and to a lesser extent with liver (HP 9314). Strong bands have also been detected in extracts of whole rat brain and HepG2, a human hepatocellular carcinoma cell line (ATCC HB8065). A single band at ~36kDa is observed at a primary antibody dilution of 1:2000 or above when detected using enhanced chemiluminescence. Antibody HC 3002 is believed to be wholly specific for HO-2 showing no cross-reactivity with HO-1 in all systems studied to date. The immunostaining is fully abolished by pre-adsorption of the antibody with cognate peptide (HP 9302). The antibody has been used successfully in immunocytochemical studies of HO-2 localisation in human bronchial epithelium biopsy material.

Fixatives: Several fixative solutions may be used. Aldehyde-combination fixatives (i.e. those containing formaldehyde and glutaraldehyde) usually give satisfactory results. Bouin and Susa fixatives containing 0.1-0.2% glutaraldehyde have been used satisfactorily.

P30519

This antibody is covered by our Worry-Free Guarantee

.

UniProt ID

Worry-free Guarantee

Last modified: May 29, 2024

ENZO LIFE SCIENCES,

ENZO LIFE SCIENCES, INC.
Phone: 800.942.0430
infousa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com