

HO1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # H20-30G

Lot # V334-3

Product Description

Recombinant full-length human HO1 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_002133](#).

Gene Aliases

HMOX1, HSP32, bK286B10

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

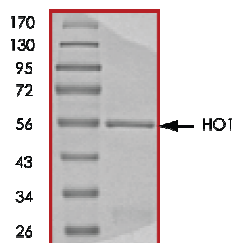
Scientific Background

HO1 (Heme oxygenase 1) catalyzes the oxidative cleavage of heme to biliverdin and is one of the main genes controlling heme synthesis and catabolism. **HO1** plays a protective role in various disorders and can ameliorate experimental MN via multiple pathways, including anti-oxidative and immunomodulatory effects (1). Exposure of primary hepatocytes to carbon monoxide and nitric oxide results in dramatic induction of HO1 in dose- and time-dependent manner and this induction is blocked by MAP kinase inhibitors (MAPKs) but not by inhibitors of PI3 kinase pathway (2).

References

1. Chia-Chao, W.u. et al: HO-1 induction ameliorates experimental murine membranous nephropathy: anti-oxidative, anti-apoptotic and immunomodulatory effects. *Nephrology Dialysis Transplantation*, 2008; 23(10):3082-3090.
2. Lee, B.S. et al: Carbon monoxide mediates heme oxygenase 1 induction via Nrf2 activation in hepatoma cells. *Biochem Biophys Res Commun*. 2006 May 12;343(3):965-72.

Purity



The purity of HO1 was determined to be **>95%** by densitometry. Approx. MW **56kDa**.

HO1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	H20-30G
Specific Lot Number	V334-3
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.