

HO2 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # H21-30G

Lot # D038-3

Product Description

Recombinant full-length human HO2 Protein was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_002134](#).

Gene Aliases

HMOX2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 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.

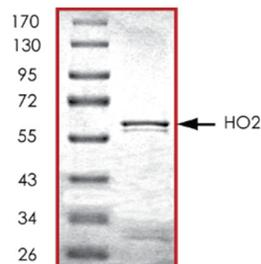
Scientific Background

HO2, or heme oxygenase 2, belongs to the heme oxygenase family that are essential enzyme in heme catabolism that cleave heme to form biliverdin which is subsequently converted to bilirubin by biliverdin reductase (1). HO2 is part of the calcium-sensitive potassium channel complex and enhances channel activity in normoxia (2). Knockdown of HO2 expression reduces channel activity, an effect that is reversed by carbon monoxide. Inhibition of calcium-sensitive potassium channels by hypoxia is dependent on HO2 expression and is augmented by HO2 stimulation.

References

1. Kemp, P.J. et al: Hemeoxygenase-2 as an O2 sensor in K+ channel-dependent chemotransduction. *Biochem Biophys Res Commun.* 2005;338(1):648-52.
2. Williams, S. E. et al: Hemoxygenase-2 is an oxygen sensor for a calcium-sensitive potassium channel. *Science.* 2004 Dec 17;306(5704):2093-7.

Purity



The purity of HO2 protein was determined to be **>85%** by densitometry. Approx. MW **62kDa**.

HO2 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog #	H21-30G
Lot #	D038-3
Purity	>85%
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.