

PRODUCT INFORMATION



HO-1 (human, recombinant)

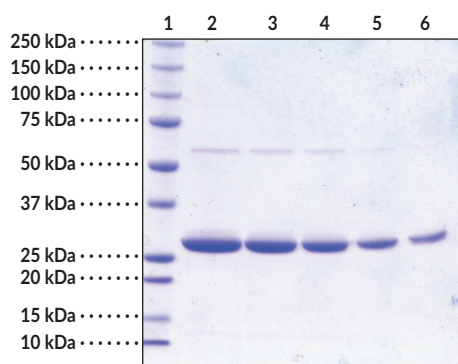
Item No. 22731

Overview and Properties

Synonyms: Heat Shock Protein 32, Heme Oxygenase-1, HMOX1, Hsp32
Source: Recombinant C-terminal His-tagged expressed in *E. coli*
Amino acids: 1-265 (full length)
Uniprot No.: P09601
Molecular Weight: 31 kDa
Storage: -80°C (as supplied)
Stability: ≥3 years
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 50 mM HEPES, pH 8.0, with 150 mM sodium chloride
Protein Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: HO-1 (8 µg)
Lane 3: HO-1 (6 µg)
Lane 4: HO-1 (4 µg)
Lane 5: HO-1 (2 µg)
Lane 6: HO-1 (1 µg)

Representative gel image shown; actual purity may vary between each batch.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/06/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Heme oxygenase-1 (HO-1), also known as heat shock protein 32 (Hsp32), is an inducible heme oxygenase encoded by the HMOX1 gene.¹⁻³ It is a membrane-bound enzyme that catalyzes the cleavage of heme to release carbon monoxide (CO), ferrous ions (Fe²⁺), and biliverdin, with biliverdin being further processed into bilirubin. HO-1 is found in human spleen, liver, and kidney where its expression is induced by the presence of heme, hormones, metals, oxidative agents, and therapeutic compounds to protect against oxidative stress and inflammatory responses. HO-1 is upregulated in a variety of cancers and siRNA knockdown of HMOX1 or inhibition of HO-1 decreases cancer cell proliferation.⁴⁻⁶ HO-1 also interacts with the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) accessory protein Orf3a that, in a similar virus, SARS-CoV, is associated with activation of the NLRP3 inflammasome.⁷⁻⁹ Cayman's HO-1 protein can be used for ELISA and Western blot (WB) applications.

References

1. Wang, J., Lad, L., Poulos, T.L., *et al.* Regiospecificity determinants of human heme oxygenase. *J. Biol. Chem.* **280(4)**, 2797-2806 (2005).
2. Martasek, P., Solangi, K., Goodman, A.I., *et al.* Properties of human kidney heme oxygenase: Inhibition by synthetic heme analogues and metalloporphyrins. *Biochem. Biophys. Res. Commun.* **157(2)**, 480-487 (1988).
3. Huber, W.J., III, Marohnic, C.C., Peters, M., *et al.* Measurement of membrane-bound human heme oxygenase-1 activity using a chemically defined assay system. *Drug Metab. Dispos.* **37(4)**, 857-864 (2009).
4. Berberat, P.O., Dambrauskas, Z., Gulbinas, A., *et al.* Inhibition of heme oxygenase-1 increases responsiveness of pancreatic cancer cells to anticancer treatment. *Clin. Cancer Res.* **11(10)**, 3790-3798 (2005).
5. Nowis, D., Legat, M., Grzela, T., *et al.* Heme oxygenase-1 protects tumor cells against photodynamic therapy-mediated cytotoxicity. *Oncogene* **25(24)**, 3365-3374 (2006).
6. Fang, J., Akaike, T., and Maeda, H. Antiapoptotic role of heme oxygenase (HO) and the potential of HO as a target in anticancer treatment. *Apoptosis* **9(1)**, 27-35 (2004).
7. Gordon, D.E., Jang, G.M., Bouhaddou, M. *et al.* A SARS-CoV-2-human protein-protein interaction map reveals drug targets and potential drug-repurposing. *BioRxiv* (2020).
8. Li, H., Liu, S.-M., Yu, X.-H. *et al.* Coronavirus disease 2019 (COVID-19): Current status and future perspectives. *Int. J. Antimicrob. Agents* (2020).
9. Siu, K.-L., Yuen, K.-S., Castaño-Rodríguez, C., *et al.* Severe acute respiratory syndrome coronavirus ORF3a protein activates the NLRP3 inflammasome by promoting TRAF3-dependent ubiquitination of ASC. *FASEB J.* **33(8)**, 8865-8877 (2019).