

PRODUCT INFORMATION



HO-2 (human recombinant)

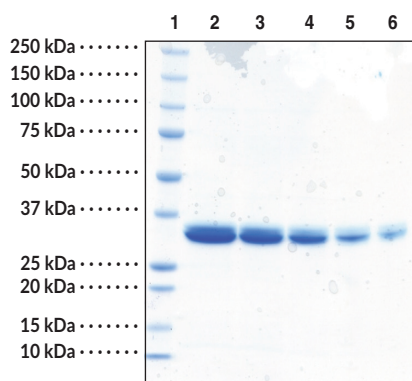
Item No. 22732

Overview and Properties

Synonyms: EC:1.14.14.18, Heme oxygenase 2, HMOX2
Source: N-terminal His-tagged HO-2/HMOX2 human recombinant protein purified from *E. coli*
Amino Acids: 2-264
Uniprot No.: P30519
Molecular Weight: 32.6 kDa
Storage: -80°C (as supplied); avoid freeze/thaw cycles by storing protein in aliquots
Stability: ≥1 year
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 50 mM HEPES, pH 8.0, 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-2 (8 µg)
Lane 3: HO-2 (6 µg)
Lane 4: HO-2 (4 µg)
Lane 5: HO-2 (2 µg)
Lane 6: HO-2 (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, 03/06/2019

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-2 (HO-2) is a constitutively active heme oxygenase encoded by the *HMOX2* gene.^{1,2} It is a membrane-bound enzyme that catalyzes the cleavage of heme to give carbon monoxide (CO), ferrous ions (Fe^{2+}), and biliverdin, with biliverdin being further processed into bilirubin. HO-2 is found in neurons, testes, and endothelial and smooth muscle cells from cerebral vessels.² HO-2 protects against apoptotic neuronal cell death in models of ischemic injury and oxidative stress. It also acts as an oxygen sensor to inhibit systemic vascular reactivity and reduce cell death in response to hypoxia.

References

1. Maines, M.D. The heme oxygenase system: A regulator of second messenger gases. *Annu. Rev. Pharmacol. Toxicol.* **37**, 517-554 (1997).
2. Muñoz-Sánchez, J. and Cháñez-Cárdenas, M.E. A review on hemeoxygenase-2: Focus on cellular protection and oxygen response. *Oxid. Med. Cell. Longev.* **604981**, (2014).

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