

PRODUCT INFORMATION



Soluble Epoxide Hydrolase (mouse, recombinant)

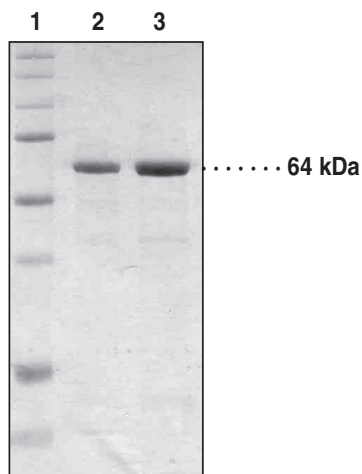
Item No. 10011670

Overview and Properties

Synonyms: Cytosolic Epoxide Hydrolase (CEH), EPHX2, Epoxide Hydrolase 2, sEH
Source: Active recombinant mouse N-terminal His-tagged sEH expressed in insect cells
Uniprot No.: P34914
Molecular Weight: 64 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: *batch specific* (≥95% estimated by SDS-PAGE)
Supplied in: TBS, pH 7.4 with 20% glycerol
Protein
Concentration: *batch specific* mg/ml
Activity: *batch specific* U/ml
Specific Activity: *batch specific* U/mg
Unit Definition: One unit is defined as the amount of enzyme required to produce 1 nmole of 6-methoxy-2-naphthaldehyde per minute at 37°C in 25 mM bis-Tris, pH 7.0, containing 1 μM PHOME

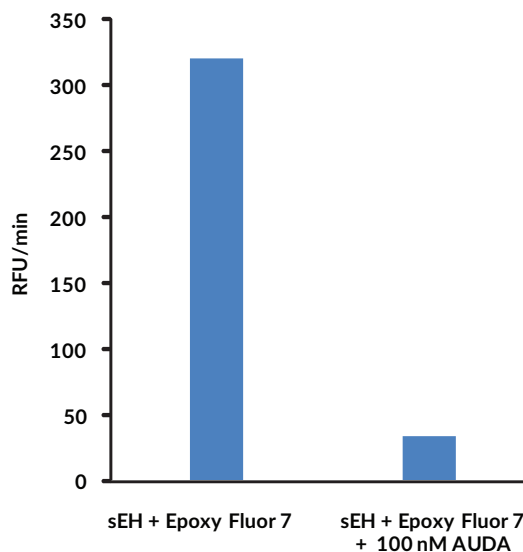
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Standards
Lane 2: Mouse sEH (2 μg)
Lane 3: Mouse sEH (4 μg)

SDS-PAGE Analysis of Soluble Epoxide Hydrolase



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, 09/14/2022

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

Soluble epoxide hydrolase (sEH) is a member of the α/β -hydrolase fold enzyme family that catalyzes the hydrolysis of bioactive fatty acid epoxides to inactive vicinal diols.¹ It is a homodimer in which each subunit is composed of two domains, a C-terminal epoxide hydrolase domain and an N-terminal phosphatase domain.^{2,3} sEH is localized to the cytoplasm or to peroxisomes in a tissue-specific manner and is found in various tissues, including skin, lung, uterus, kidney, brain, and myocardium.^{4,5} sEH is also expressed in the vasculature and inhibition of sEH attenuates pathogenic vascular remodeling and hypertension via preservation of cardioprotective epoxyeicosatrienoic acids (EETs) in rat models of atherosclerosis and hypertension, respectively.⁶ Inhibition of sEH also has a protective role in various diseases, including inflammatory bowel disease, osteoarthritis, seizure, stroke, and Alzheimer's disease, as well as in various chronic pain states.^{1,7} Cayman's Soluble Epoxide Hydrolase (mouse, recombinant) protein can be used for enzyme activity assays.

References

1. Harris, T.R. and Hammock, B.D. Soluble epoxide hydrolase: Gene structure, expression and deletion. *Gene* **526(2)**, 61-74 (2013).
2. Cronin, A., Mowbray, S., Dürk, H., et al. The N-terminal domain of mammalian soluble epoxide hydrolase is a phosphatase. *Proc. Natl. Acad. Sci. USA* **100(4)**, 1552-1557 (2002).
3. Chiamvimonvat, N., Ho, C.M., Tsai, H.J., et al. The soluble epoxide hydrolase as a pharmaceutical target for hypertension. *J. Cardiovasc. Pharmacol.* **50(3)**, 225-237 (2007).
4. Enayetallah, A.E., French, R.A., Barber, M., et al. Cell-specific subcellular localization of soluble epoxide hydrolase in human tissues. *J. Histochem. Cytochem.* **54(3)**, 329-335 (2006).
5. Domingues, M.F., Callai-Silva, N., Piovesan, A.R., et al. Soluble epoxide hydrolase and brain cholesterol metabolism. *Front. Mol. Neurosci.* **12**, 325 (2020).
6. Simpkins, A.N., Rudic, R.D., Roy, S., et al. Soluble epoxide hydrolase inhibition modulates vascular remodeling. *Am. J. Physiol. Heart Circ. Physiol.* **298(3)**, H795-H806 (2009).
7. Wagner, K.M., McReynolds, C.B., Schmidt, W.K., et al. Soluble epoxide hydrolase as a therapeutic target for pain, inflammatory and neurodegenerative diseases. *Pharmacol. Ther.* **180**, 62-76 (2017).

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