

PRODUCT INFORMATION



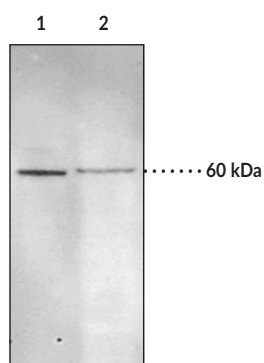
Soluble Epoxide Hydrolase Polyclonal Antibody

Item No. 10010146

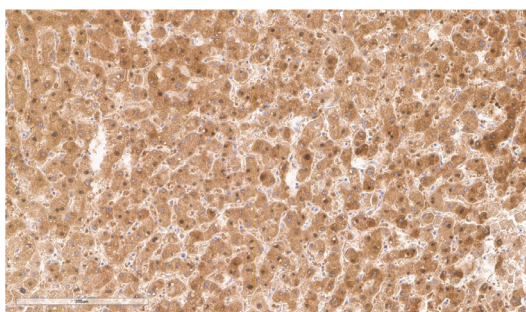
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	Cytosolic Epoxide Hydrolase (CEH), EPHX2, Epoxide Hydrolase 2, sEH
Immunogen:	Synthetic peptide from the internal region of human sEM
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	P34913
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Human liver 10,000 x g supernatant (30 µg)
Lane 2: Mouse liver 10,000 x g supernatant (60 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human liver tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with Soluble Epoxide Hydrolase Polyclonal Antibody (Item No. 10010146) at a 1:40 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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, 10/17/2023

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 Polyclonal Antibody can be used for Western blot. The antibody recognizes sEH at approximately 64 kDa from human, mouse, and rat samples.

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