

# PRODUCT INFORMATION



## Soluble Epoxide Hydrolase Blocking Peptide

Item No. 10010147

### Overview and Properties

---

<b>Contents:</b>	This vial contains 200 µg of lyophilized peptide.
<b>Synonyms:</b>	Cytosolic Epoxide Hydrolase (CEH), EPHX2, Epoxide Hydrolase, sEH
<b>Form:</b>	Solid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years

### Procedures

---

This vial contains 200 µg of lyophilized peptide. This blocking peptide can be used in conjunction with soluble epoxide hydrolase polyclonal antibody (Item No. 10010146) to block protein-antibody complex formation during immunochemical analysis of soluble epoxide hydrolase.

To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the Soluble Epoxide Hydrolase Polyclonal Antibody (Item No. 10010146) at the recommended concentration (1:200 for Western blot (WB)) with 10 µg/ml blocking peptide at the volume desired for the application (10 ml typical for WB).\*
2. Incubate for one hour at room temperature with occasional mixing.
3. After one hour of incubation, apply to the slide or membrane as usual.

\*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

#### 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, 04/04/2024

#### 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