

Product datasheet

Soluble Epoxide hydrolase Inhibitor Screening Kit (Fluorometric) ab283390

[2 Images](#)

Overview

Product name	Soluble Epoxide hydrolase Inhibitor Screening Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Inhibitor compounds
Assay duration	Multiple steps standard assay
Product overview	Soluble Epoxide hydrolase (sEH) Inhibitor Screening Kit (Fluorometric) (ab283390) can be used to screen/characterize/study potential inhibitors of Soluble Epoxide hydrolase.

This kit is a micro-plate based fluorometric kit for screening sEH inhibitors. Its assay principle is based on the hydrolysis of a non-fluorescent substrate to a fluorescent product by sEH. It uses N-Cyclohexyl-N'-dodecylurea (NCND), a well-known sEH inhibitor, as positive control.

The assay is rapid, simple and sensitive.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K480 Soluble Epoxide Hydrolase Inhibitor Screening Kit (Fluorometric). K480-100 is the same size as the 100 test size of ab283390.

Platform Microplate (12 x 8 well strips)

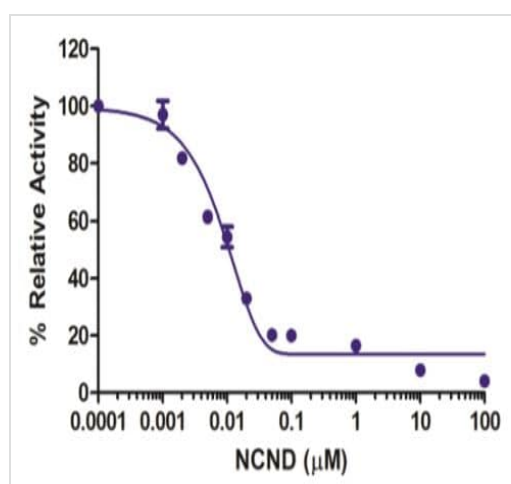
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Epoxide hydrolase Assay Buffer	1 x 25ml
Epoxide hydrolase Substrate	1 x 200µl
Human Epoxide hydrolase Enzyme	1 vial
NCND (5 mM)	1 x 100µl

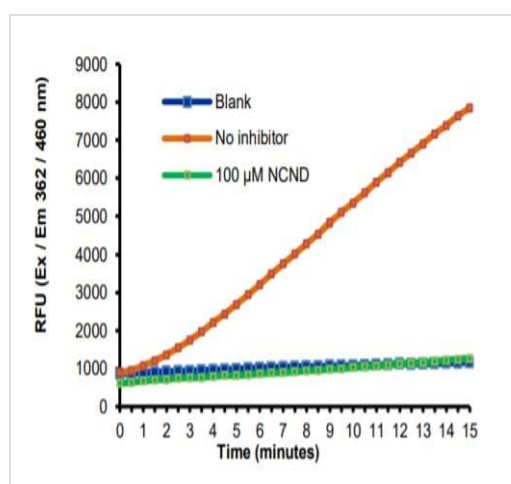
Function	Biotransformation enzyme that catalyzes the hydrolysis of arene and aliphatic epoxides to less reactive and more water soluble dihydrodiols by the trans addition of water.
Tissue specificity	Found in liver.
Involvement in disease	Note=In some populations, the high activity haplotype tyr113/his139 is overrepresented among women suffering from pregnancy-induced hypertension (pre-eclampsia) when compared with healthy controls. Defects in EPHX1 are a cause of familial hypercholanemia (FHCA) [MIM:607748]. FHCA is a disorder characterized by elevated serum bile acid concentrations, itching, and fat malabsorption.
Sequence similarities	Belongs to the peptidase S33 family.
Cellular localization	Microsome membrane. Endoplasmic reticulum membrane.

Images



Inhibition of soluble Epoxide Hydrolase enzyme activity by NCND.
IC50 of NCND was determined to be 7.15 ± 0.43 nM.

Functional Studies - Soluble Epoxide hydrolase
Inhibitor Screening Kit (Fluorometric) (ab283390)



sEH enzyme kinetics in presence and absence of Inhibitor NCND.

Functional Studies - Soluble Epoxide hydrolase
Inhibitor Screening Kit (Fluorometric) (ab283390)

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors