

# PRODUCT INFORMATION



## Soluble Epoxide Hydrolase PROTAC 1a

Item No. 38876

**Formal Name:** N-(1-((2-(2,6-dioxopiperidin-3-yl)-1,3-dioxisoindolin-4-yl)oxy)-2-oxo-6,9,12,15-tetraoxa-3-azaheptadecan-17-yl)-4-(((1r,4r)-4-(3-(4-(trifluoromethoxy)phenyl)ureido)cyclohexyl)oxy)benzamide

**Synonyms:** sEH PROTAC 1a, sEH Proteolysis-targeting Chimera 1a, Soluble Epoxide Hydrolase Proteolysis-targeting Chimera 1a

**MF:** C<sub>46</sub>H<sub>53</sub>F<sub>3</sub>N<sub>6</sub>O<sub>14</sub>

**FW:** 971.0

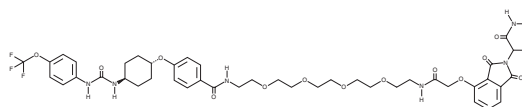
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 245 nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥3 years



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

### Description

Soluble epoxide hydrolase (sEH) PROTAC 1a is a proteolysis-targeting chimera (PROTAC) that contains cereblon ligand 1 (Item No. 21589) connected to the sEH inhibitor *t*-TUCB (Item No. 36384) by a linker region.<sup>1</sup> It drives sEH degradation and selectively inhibits the hydrolase activity of sEH (IC<sub>50</sub> = 0.8 nM) over its phosphatase activity (IC<sub>50</sub> = >10,000 nM). sEH PROTAC 1a is also selective for the degradation of cytosolic over peroxisomal sEH and drives its degradation *via* the lysosome rather than the proteasome. It reduces thapsigargin-induced increases in phosphorylated inositol-requiring enzyme 1α (IRE1α) levels and X-box binding protein 1 (XBP1) splicing in HepG2 and 293T cells, indicating a reduction in ER stress.

### Reference

1. Wang, Y., Morisseau, C., Takamura, A., *et al.* PROTAC-Mediated selective degradation of cytosolic soluble epoxide hydrolase enhances ER stress reduction. *ACS Chem. Biol.* **18**(4), 884-896 (2023).

#### 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, 07/12/2023

#### 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