

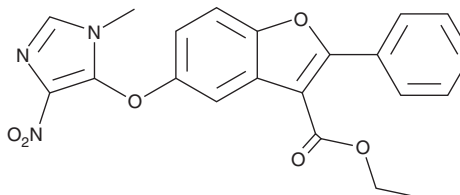
# PRODUCT INFORMATION



## STOML3 Inhibitor OB-1

Item No. 37584

**CAS Registry No.:** 300803-69-4  
**Formal Name:** 5-[(1-methyl-4-nitro-1H-imidazol-5-yl)oxy]-2-phenyl-3-benzofurancarboxylic acid, ethyl ester  
**Synonym:** Stomatin-like Protein 3 Inhibitor OB-1  
**MF:** C<sub>21</sub>H<sub>17</sub>N<sub>3</sub>O<sub>6</sub>  
**FW:** 407.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 306 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

STOML3 inhibitor OB-1 is supplied as a solid. A stock solution may be made by dissolving the STOML3 inhibitor OB-1 in the solvent of choice, which should be purged with an inert gas. STOML3 inhibitor OB-1 is soluble in DMSO.

### Description

STOML3 inhibitor OB-1 is an inhibitor of stomatin-like protein 3 (STOML3) oligomerization.<sup>1</sup> It inhibits oligomerization of mouse and human STOML3 in cell-free assays when used at a concentration of 10 μM. STOML3 inhibitor OB-1 reduces Piezo1 channel current amplitude in Neuro2a cells. *In vivo*, STOML3 inhibitor OB-1 inhibits tactile perception in wild-type mice and reduces tactile allodynia in a mouse model of diabetic neuropathy induced by streptozotocin (Item No. 13104).

### Reference

1. Wetzel, C., Pifferi, S., Picci, C., *et al.* Small-molecule inhibition of STOML3 oligomerization reverses pathological mechanical hypersensitivity. *Nat. Neurosci.* **20(2)**, 209-218 (2017).

#### 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, 11/14/2022

#### 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