

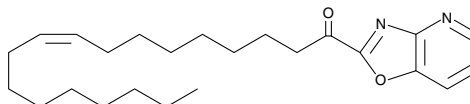
# Product Information



## Oleoyl Oxazolopyridine

Item No. 71650

**CAS Registry No.:** 288862-58-8  
**Formal Name:** 1-oxazolo[4,5-b]pyridin-2-yl-octadec-9Z-en-1-one  
**Synonym:** CAY10400  
**MF:** C<sub>24</sub>H<sub>36</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 384.6  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in methyl acetate  
**UV/Vis.:** λ<sub>max</sub>: 234, 297 nm



### Laboratory Procedures

For long term storage, we suggest that oleoyl oxazolopyridine be stored as supplied at -20°C. It should be stable for at least one year.

Oleoyl oxazolopyridine is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of oleoyl oxazolopyridine in these solvents is approximately 10 mg/ml.

Oleoyl oxazolopyridine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, first evaporate the methyl acetate under a gentle stream of nitrogen and dissolve the neat oil in ethanol. The ethanolic solution can then be diluted with the aqueous buffer of choice. Oleoyl oxazolopyridine has a solubility of approximately 50 µg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Fatty acid amide hydrolase (FAAH) is the enzyme responsible for hydrolysis and inactivation of fatty acid amides including anandamide and oleamide. CAY10400 is a potent inhibitor of FAAH exhibiting K<sub>i</sub> values of 1.3 nM and 2.3 nM for the human and rat enzymes, respectively.<sup>1</sup> Oleoyl oxazolopyridine is approximately 35-fold more potent than oleyl trifluoromethyl ketone when assayed under the same conditions. Selectivity of CAY10400 for FAAH *versus* cannabinoid receptors or PLA<sub>2</sub> has not been reported.

### Reference

1. Boger, D.L., Sato, H., Lerner, A.E., *et al.* Exceptionally potent inhibitors of fatty acid amide hydrolase: The enzyme responsible for degradation of endogenous oleamide and anandamide. *Proc. Natl. Acad. Sci. USA* **97**, 5044-5049 (2000).

### Related Products

URB597 - Item No. 10046 • N-Oleoyl Dopamine - Item No. 10115 • URB937 - Item No. 10674 • Oleyl Trifluoromethyl Ketone - Item No. 62640 • Arachidonoyl Serotonin - Item No. 70665

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/09/2011

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com