

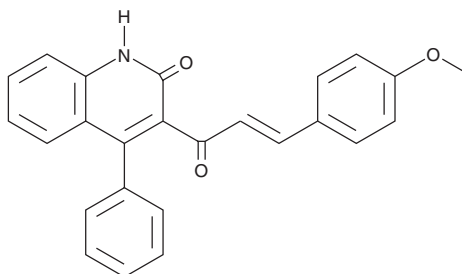
PRODUCT INFORMATION



Ceranib-2

Item No. 11092

CAS Registry No.: 1402830-75-4
Formal Name: 3-[3-(4-methoxyphenyl)-1-oxo-2-propen-1-yl]-4-phenyl-2(1H)-quinolinone
MF: C₂₅H₁₉NO₃
FW: 381.4
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 338 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Ceranib-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the ceranib-2 in the solvent of choice. Ceranib-2 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ceranib-2 in DMSO is approximately 20 mg/ml and approximately 25 mg/ml in DMF.

Ceranib-2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ceranib-2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ceranib-2 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ceranib-2 is a ceramidase inhibitor that inhibits ceramidase activity (IC₅₀ = 28 μM in SKOV3 cells).¹ It also increases intracellular accumulation of various ceramides, including C14 ceramide (Item No. 22531), C24 ceramide (Item No. 62535), C26 ceramide, and C26:1 ceramide, and total ceramide levels and decreases levels of dihydrosphingosine, sphingosine, and sphingosine-1-phosphate (Item No. 22500) in SKOV3 cells. *In vivo*, ceranib-2 (20 and 50 mg/kg, i.p.) reduces tumor volume in a JC mammary adenocarcinoma mouse syngeneic tumor model.

Reference

1. Draper, J.M., Xia, Z., Smith, R.A., *et al.* Discovery and evaluation of inhibitors of human ceramidase. *Mol. Cancer Ther.* **1-27** (2011).

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, 07/19/2018

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