

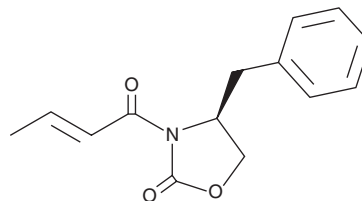
PRODUCT INFORMATION



(E/Z)-Locostatin

Item No. 29092

CAS Registry No.: 133812-16-5
Formal Name: (4S)-3-(1-oxo-2-buten-1-yl)-4-(phenylmethyl)-2-oxazolidinone
MF: $C_{14}H_{15}NO_3$
FW: 245.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 213, 217, 226 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

(E/Z)-Locostatin is supplied as a crystalline solid. A stock solution may be made by dissolving the (E/Z)-locostatin in the solvent of choice, which should be purged with an inert gas. (E/Z)-Locostatin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (E/Z)-locostatin in ethanol is approximately 10 mg/ml and approximately 5 mg/ml in DMSO and DMF.

(E/Z)-Locostatin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (E/Z)-locostatin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (E/Z)-Locostatin has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(E/Z)-Locostatin is the racemic form of locostatin, an inhibitor of cell migration and the protein-protein interaction between Raf kinase inhibitor protein (RKIP) and Raf-1.^{1,2}

References

1. Mc Henry, K.T., Ankala, S.V., Ghosh, A.K., *et al.* A non-antibacterial oxazolidinone derivative that inhibits epithelial cell sheet migration. *Chembiochem.* **3(11)**, 1105-1111 (2002).
2. Zhu, S., Mc Henry, K.T., Lane, W.S., *et al.* A chemical inhibitor reveals the role of Raf kinase inhibitor protein in cell migration. *Chem. Biol.* **12(9)**, 981-991 (2005).

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, 11/16/2022

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