

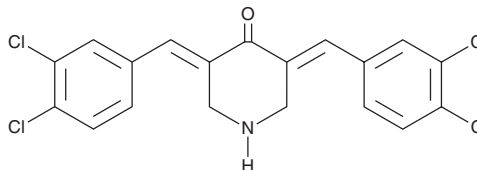
PRODUCT INFORMATION



RAMB4

Item No. 36936

CAS Registry No.: 145888-79-5
Formal Name: 3,5-bis[(3,4-dichlorophenyl)methylene]-4-piperidinone
Synonym: NSC 638643
MF: C₁₉H₁₃Cl₄NO
FW: 413.1
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 329 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

RAMB4 is supplied as a solid. A stock solution may be made by dissolving the RAMB4 in the solvent of choice, which should be purged with an inert gas. RAMB4 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

RAMB4 is an inhibitor of the ubiquitin-proteasome system.¹ It increases the levels of polyubiquitinated proteins in HeLa cervical cancer cells, but does not inhibit the chymotrypsin-like, trypsin-like, or peptidylglutamyl peptide hydrolyzing-like activities of the 20S proteasome, in cell-free assays when used at a concentration of 10 μM. RAMB4 decreases the viability of HeLa cells (IC₅₀ = 2 μM).

Reference

1. Anchoori, R.K., Khan, S.R., Sueblinvong, T., *et al.* Stressing the ubiquitin-proteasome system without 20S proteolytic inhibition selectively kills cervical cancer cells. *PLoS One* **6**(8), e23888 (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, 10/31/2022

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