

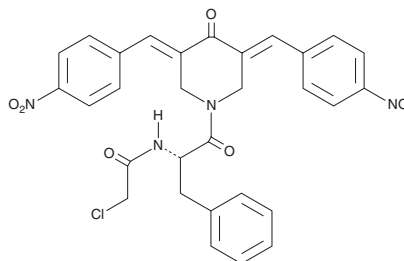
PRODUCT INFORMATION



RA-375

Item No. 41541

CAS Registry No.: 2649154-57-2
Formal Name: N-[(1S)-2-[(3E,5E)-3,5-bis[(4-nitrophenyl)methylene]-4-oxo-1-piperidiny]-2-oxo-1-(phenylmethyl)ethyl]-2-chloro-acetamide
MF: C₃₀H₂₅ClN₄O₇
FW: 589.0
Purity: ≥98%
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

RA-375 is supplied as a solid. A stock solution may be made by dissolving the RA-375 in the solvent of choice, which should be purged with an inert gas. RA-375 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

RA-375 is an inhibitor of the ubiquitin-proteasome system.¹ It inhibits the degradation of a polyubiquitinated reporter protein in a reporter assay using HEK293T cells when used at concentrations of 0.3125 or 0.625 μ M. RA-375 (1 μ M) increases levels of mRNA encoding DNA damage-inducible transcript 3 (DDIT3), also known as C/EBP homologous protein (CHOP), and X-box binding protein 1 (XBP1), both of which are endoplasmic reticulum (ER) stress markers, in ES-2 ovarian cancer cells. It also increases levels of reactive oxygen species (ROS) and induces apoptosis in ES-2 cells when used at a concentration of 1 μ M. *In vivo*, RA-375 reduces tumor growth in an ES-2 mouse xenograft model.

Reference

1. Anchoori, R.K., Tan, M., Tseng, S.H., *et al.* Structure-function analyses of candidate small molecule RPN13 inhibitors with antitumor properties. *PLoS One* **15**(1), e0227727 (2020).

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, 09/18/2024

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