

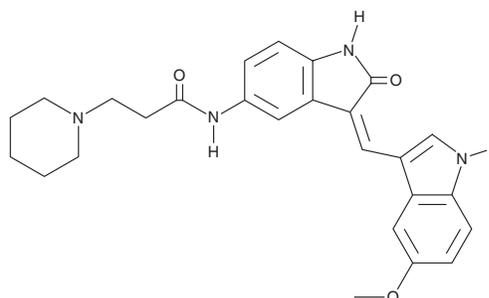
PRODUCT INFORMATION



DEL-22379

Item No. 22433

CAS Registry No.: 181223-80-3
Formal Name: N-[2,3-dihydro-3-[(5-methoxy-1H-indol-3-yl)methylene]-2-oxo-1H-indol-5-yl]-1-piperidinepropanamide
MF: C₂₆H₂₈N₄O₃
FW: 444.5
Purity: ≥98% (mixture of rotamers)
UV/Vis.: λ_{max}: 224, 276, 405 nm
Supplied as: A crystalline 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

DEL-22379 is supplied as a crystalline solid. A stock solution may be made by dissolving the DEL-22379 in the solvent of choice, which should be purged with an inert gas. DEL-22379 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of DEL-22379 in these solvents is approximately 30 mg/ml. DEL-22379 is also slightly soluble in ethanol.

DEL-22379 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DEL-22379 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DEL-22379 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DEL-22379 is an inhibitor of ERK dimerization that disrupts EGF-induced co-immunoprecipitation of ERK tagged with hemagglutinin (HA) or FLAG epitopes with an IC₅₀ value of approximately 0.5 μM without affecting ERK phosphorylation or kinase activity (IC₅₀s = >10 μM).¹ It inhibits cell growth in a panel of melanoma and colorectal cancer cell lines (IC₅₀s = 150-400 nM). DEL-22379 (15 mg/kg, i.p.) inhibits ERK dimerization in liver extract and tumor tissue and inhibits tumor progression in an A375 mouse xenograft model. It also reduces tumor growth in a BRAF mutant colorectal cancer patient-derived xenograft (PDX) mouse model.

Reference

1. Herrero, A., Pinto, A., Colón-Bolea, P., *et al.* Small molecule inhibition of ERK dimerization prevents tumorigenesis by RAS-ERK pathway oncogenes. *Cancer Cell* **28(2)**, 170-182 (2015).

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