

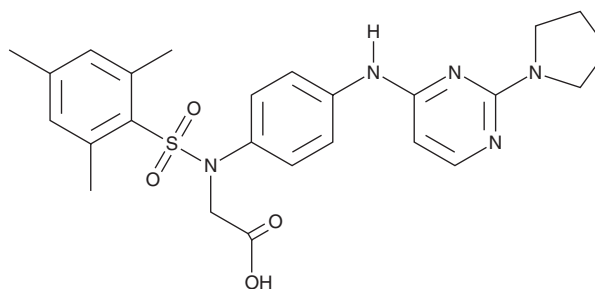
PRODUCT INFORMATION



DDO-5936

Item No. 41040

CAS Registry No.: 2355377-13-6
Formal Name: N-[4-[[2-(1-pyrrolidinyl)-4-pyrimidinyl] amino]phenyl]-N-[(2,4,6-trimethylphenyl) sulfonyl]-glycine
MF: C₂₅H₂₉N₅O₄S
FW: 495.6
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

DDO-5936 is supplied as a solid. A stock solution may be made by dissolving the DDO-5936 in the solvent of choice, which should be purged with an inert gas. DDO-5936 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in chloroform and acetonitrile.

Description

DDO-5936 is an inhibitor of the protein-protein interaction between heat shock protein 90 (Hsp90) and cell division cycle 37 (Cdc37).¹ It inhibits the Hsp90-Cdc37 interaction in HCT116 human colorectal cancer cells when used at concentrations of 5, 10, and 25 μM. DDO-5936 is selective for inhibition of the Hsp90-Cdc37 interaction over inhibition of 20 kinases (IC₅₀s = >100 μM for all) and inhibits the proliferation of HCT116 cells (IC₅₀ = 8.99 μM). It also reduces the levels of the Hsp90 kinase client Cdk4 and induces cell cycle arrest at the G₀/G₁ phase in a Cdc37-dependent manner in HCT116 cells when used at concentrations of 10 and 25 μM. DDO-5936 (50 and 100 mg/kg per day) reduces tumor growth and tumor weight in an HCT116 mouse xenograft model.

Reference

1. Wang, L., Zhang, L., Li, L., *et al.* Small-molecule inhibitor targeting the Hsp90-Cdc37 protein-protein interaction in colorectal cancer. *Sci. Adv.* **5(9)**, eaax2277 (2019).

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, 07/08/2024

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