

PRODUCT INFORMATION



GDC-6036

Item No. 41852

CAS Registry No.: 2417987-45-0

Formal Name: 1-[[[(3S)-4-[(7R)-7-[6-amino-4-methyl-3-(trifluoromethyl)-2-pyridinyl]-6-chloro-8-fluoro-2-[[[(2S)-1-methyl-2-pyrrolidinyl]methoxy]-4-quinazolinyl]-3-methyl-1-piperazinyl]-2-propen-1-one

Synonyms: Divarasib, RG-6330

MF: C₂₉H₃₂ClF₄N₇O₂

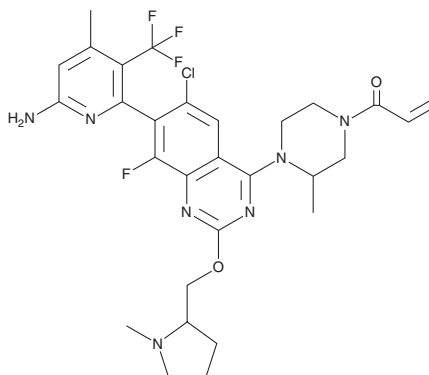
FW: 622.1

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

GDC-6036 is supplied as a solid. A stock solution may be made by dissolving the GDC-6036 in the solvent of choice, which should be purged with an inert gas. GDC-6036 is soluble (≥10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol

Description

GDC-6036 is a covalent inhibitor of K-Ras containing a glycine-to-cysteine substitution at position 12 (K-Ras^{G12C}).¹ It decreases tumor volume in a PaCa-2 pancreatic cancer mouse xenograft model when administered at doses ranging from 1 to 100 mg/kg per day.

Reference

1. Tran, J.C., Hunsaker, T., Bell, C., *et al.* Quantifying KRAS G12C covalent drug inhibitor activity in mouse tumors using mass spectrometry. *Anal. Chem.* **95**(11), 4834-4839 (2023).

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, 10/07/2024

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