

PRODUCT INFORMATION

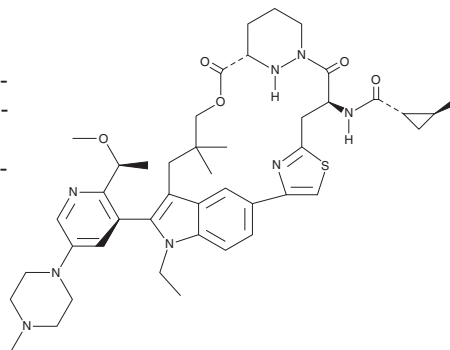


RMC-6236

Item No. 41982

CAS Registry No.: 2765081-21-6
Formal Name: (1S,2S)-N-[(2R,14S,18S)-1-ethyl-18,19,20,21-tetrahydro-2-[2-[(1S)-1-methoxyethyl]-5-(4-methyl-1-piperazinyl)-3-pyridinyl]-25,25-dimethyl-15,22-dioxo-17H-5,3-[4,2]-endo-thiazolopropano[1,3]-endo-pyridazinomethanoxypropano]-1H-indol-14-yl]-2-methyl-cyclopropanecarboxamide

MF: C₄₄H₅₈N₈O₅S
FW: 811.1
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

RMC-6236 is supplied as a solid. A stock solution may be made by dissolving the RMC-6236 in the solvent of choice, which should be purged with an inert gas. RMC-6236 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

RMC-6236 is a non-covalent inhibitor of GTP-bound RAS.¹ It binds to peptidyl-prolyl *cis-trans* isomerase A (PPIase A), also known as cyclophilin A (CypA; $K_d = 55.3$ nM) to form a binary complex that then binds to wild-type K-Ras, K-Ras(G12V), and K-Ras(G12D) ($K_{d,s} = 131, 364, \text{ and } 154$ nM, respectively). RMC-6236 forms a trimeric complex with PPIase A and a K-Ras, H-Ras, N-Ras, or mutant Ras variant and disrupts the protein-protein interaction between these Ras variants and the B-RAF Ras-binding domain ($IC_{50s} = 28-220$ nM). It induces apoptosis in 28 K-Ras(G12X) and five N-Ras(Q61X) mutant cancer cell lines when used at concentrations ranging from 1 to 100 nM. *In vivo*, RMC-6236 (25 mg/kg) inhibits intratumoral ERK phosphorylation and reduces tumor volume in Capan-2 and NCI H441 pancreatic ductal adenocarcinoma (PDAC) mouse xenograft models. It also reduces tumor growth in 29 non-small cell lung cancer (NSCLC), 22 PDAC, and 23 colorectal cancer K-Ras(G12X) mutant mouse xenograft models when administered at a dose of 25 mg/kg.

Reference

- Jiang, J., Jiang, L., Maldonato, B.J., *et al.* Translational and therapeutic evaluation of RAS-GTP inhibition by RMC-6236 in RAS-driven cancers. *Cancer Discov.* **14**(6), 994-1017 (2024).

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, 09/04/2024

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