

PRODUCT INFORMATION



RMC-6291

Item No. 42174

CAS Registry No.: 2641998-63-0

Formal Name: (3S)-N1-[N-[[1-[4-(dimethylamino)-4-methyl-1-oxo-2-pentyn-1-yl]-4-fluoro-4-piperidiny] carbonyl]-N-methyl-L-valyl-3-[4-[(2R)-1-ethyl-3-(3-hydroxy-2,2-dimethylpropyl)-2-[2-[(1S)-1-methoxyethyl]-3-pyridinyl]-1H-indol-5-yl]-2-morpholinyl]-L-alanyl]hexahydro-3-pyridazinecarboxylic acid, (3→2)-lactone

MF: C₅₅H₇₈FN₉O₈

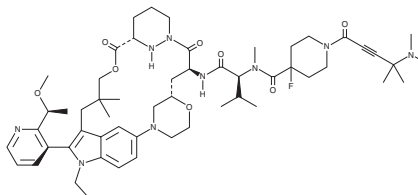
FW: 1,012.3

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-6291 is supplied as a solid. A stock solution may be made by dissolving the RMC-6291 in the solvent of choice, which should be purged with an inert gas. RMC-6291 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

RMC-6291 is a covalent inhibitor of KRAS containing a glycine-to-cysteine mutation at position 12 (KRAS^{G12C}).¹ It binds to KRAS^{G12C} and prevents KRAS^{G12C}-induced signaling via recruitment of peptidyl-prolyl *cis-trans* isomerase A (PPIase A), also known as cyclophilin A (CypA). It decreases the proliferation of 15 cancer cell lines expressing KRAS^{G12C} (median EC₅₀ = 0.11 nM) and induces apoptosis in NCI H358 non-small cell lung cancer (NSCLC) cells. RMC-6291 (10-200 mg/kg per day) reduces tumor volume in an NCI H358 mouse xenograft model.

Reference

1. Schulze, C.J., Seamon, K.J., Zhao, Y., *et al.* Chemical remodeling of a cellular chaperone to target the active state of mutant KRAS. *Science* **381**(6659), 794-799 (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, 11/12/2024

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