

PRODUCT INFORMATION



PROTAC RAF Degradar 1

Item No. 40967

CAS Registry No.: 2413035-41-1

Formal Name: (2S,4R)-1-((S)-2-(2-(4-(4-(3-(2,6-difluoro-3-(propylsulfonamido)benzoyl)-1H-pyrrolo[2,3-b]pyridin-5-yl)phenyl)piperazin-1-yl)acetamido)-3,3-dimethylbutanoyl)-4-hydroxy-N-(4-(4-methylthiazol-5-yl)benzyl)pyrrolidine-2-carboxamide

Synonyms: Proteolysis-targeting Chimera RAF Degradar 1, SJF-0628

MF: $C_{51}H_{57}F_2N_9O_7S_2$

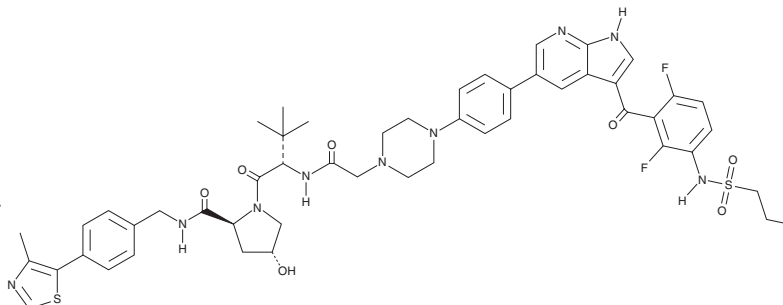
FW: 1,010.2

Purity: $\geq 98\%$

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

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

Description

PROTAC RAF degrader 1 is a proteolysis-targeting chimera (PROTAC) that contains the inhibitor of mutant V600E and wild-type B-RAF vemurafenib (PLX4032; Item No. 10618) conjugated to VHL ligand 1 (Item No. 21591).¹ It selectively induces the degradation of B-RAF containing a valine-to-glutamate substitution at position 600 (B-RAF^{V600E}) over wild-type B-RAF in SK-MEL-239 melanoma cells at 10 to 1,000 nM but non-selectively inhibits the activity of wild-type B-RAF and a panel of eight mutant B-RAF isoforms (IC_{50} s = 5.8 and 1.87-16 nM). PROTAC RAF degrader 1 reduces cell viability of SK-MEL-28, SK-MEL-239, and SK-MEL-246 melanoma cells (EC_{50} s = 215, 37, and 243 nM, respectively). *In vivo*, PROTAC RAF degrader 1 (50 mg/kg twice per day) reduces tumor growth and volume in an SK-MEL-246 mouse xenograft model.

Reference

1. Alabi, S., Jaime-Figueroa, S., Yao, Z., *et al.* Mutant-selective degradation by BRAF-targeting PROTACs. *Nat. Commun.* **12**(1), 920 (2021).

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, 03/06/2026

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