

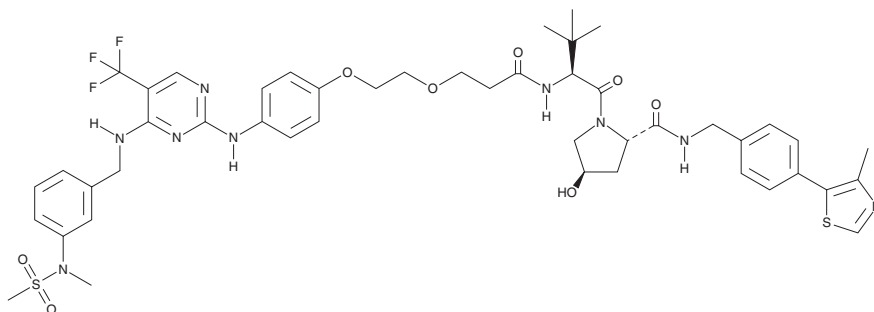
PRODUCT INFORMATION



PROTAC FAK Degradar-1

Item No. 45149

CAS Registry No.: 2301916-69-6
Formal Name: (4R)-3-methyl-N-[3-[2-[4-[[4-[[[3-[methyl(methylsulfonyl)amino]phenyl)methyl]amino]-5-(trifluoromethyl)-2-pyrimidinyl]amino]phenoxy]ethoxy]-1-oxopropyl]-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide
Synonyms: PROTAC Focal Adhesion Kinase Degradar-1 Proteolysis-targeting Chimera FAK Degradar-1 Proteolysis-targeting Chimera Focal Adhesion Kinase Degradar-1
MF: C₄₇H₅₆F₃N₉O₈S₂
FW: 996.1
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Description

PROTAC FAK degrader-1 is a proteolysis-targeting chimera (PROTAC) containing a derivative of the focal adhesion kinase (FAK) inhibitor defactinib (Item No. 17737) conjugated to VHL ligand 1 (Item No. 21591) via an alkyl ether linker.¹ PROTAC FAK degrader-1 reduces the migration and invasion of MDA-MB-231 cells when used at concentrations of 50 and 100 nM, respectively.

Reference

1. Cromm, P. M., Samarasinghe, K. T. G., Hines, J., *et al.* Addressing kinase-independent functions of fak via PROTAC-mediated degradation. *J. Am. Chem. Soc.* **140(49)**, 17019-17026 (2018).

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, 05/04/2026

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