

PRODUCT INFORMATION



Homer

Item No. 41325

CAS Registry No.: 2737222-94-3

Formal Name: N-(4'-((5-(((S)-1-((2S,4R)-4-hydroxy-2-((4-(4-methylthiazol-5-yl)benzyl)carbamoyl)pyrrolidin-1-yl)-3,3-dimethyl-1-oxobutan-2-yl)amino)-5-oxopentyl)carbamoyl)-4-(4-methylpiperazin-1-yl)-[1,1'-biphenyl]-3-yl)-6-oxo-4-(trifluoromethyl)-1,6-dihydropyridine-3-carboxamide

MF: C₅₂H₆₀F₃N₉O₇S

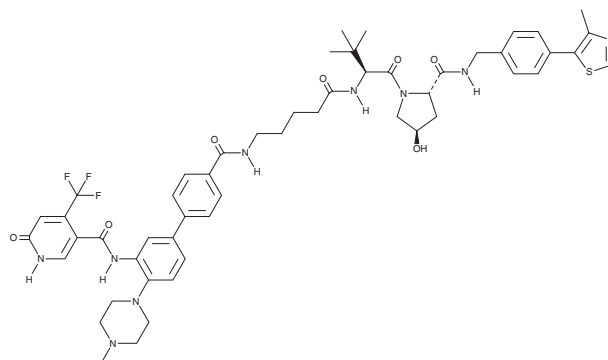
FW: 1,012.2

Purity: ≥95%

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

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

Description

Homer is a proteolysis-targeting chimera (PROTAC) composed of the probe for WD repeat domain 5 (WDR5) OICR-9429 (Item No. 16095) and a von Hippel-Lindau (VHL) ligand.¹ It induces degradation of WDR5 in MV4-11 cells with a half-maximal degradation concentration (DC₅₀) of 53 nM. Homer (1 μM) reduces WDR5 protein levels, but not levels of mRNA encoding WDR5, in MV4-11 cells. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Dölle, A., Adhikari, B., Krämer, A., *et al.* Design, synthesis, and evaluation of WD-Repeat-Containing Protein 5 (WDR5) degraders. *J. Med. Chem.* **64**(15), 10682-10710 (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, 08/09/2024

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