

PRODUCT INFORMATION



ERD-308

Item No. 33388

CAS Registry No.: 2320561-35-9

Formal Name: (2S,4R)-1-((S)-2-(2-((5-(ethyl(2-(4-(6-hydroxy-2-(4-hydroxyphenyl)benzo[b]thiophene-3-carbonyl)phenoxy)ethylamino)pentyl)oxy)acetamido)-3,3-dimethylbutanoyl)-4-hydroxy-N-((S)-1-(4-(4-methylthiazol-5-yl)phenyl)ethyl)pyrrolidine-2-carboxamide

MF: C₅₅H₆₅N₅O₉S₂

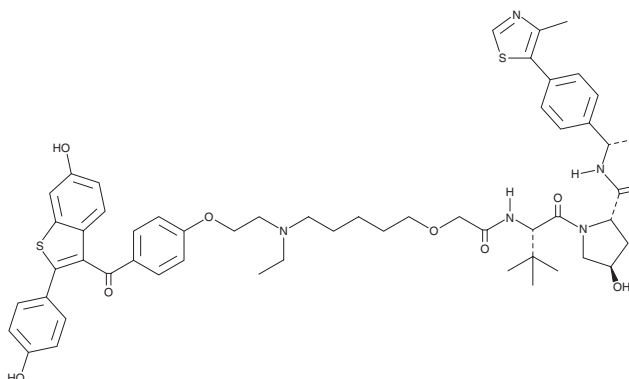
FW: 1,004.3

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

ERD-308 is supplied as a solid. A stock solution may be made by dissolving the ERD-308 in the solvent of choice, which should be purged with an inert gas. ERD-308 is slightly soluble (0.1-1 mg/ml) in methanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

ERD-308 is a proteolysis-targeting chimera (PROTAC) containing the selective estrogen receptor modulator raloxifene (Item No. 10011620) conjugated to VHL ligand 1 (Item No. 21591).¹ It induces degradation of estrogen receptor α (ER α) in MCF-7 and T47D breast cancer cells with half-maximal degradation concentration (DC₅₀) values of 0.17 and 0.43 nM, respectively. ERD-308 inhibits the growth of MCF-7 cells (GI₅₀ = 0.77 nM). It reduces the levels of mRNA encoding progesterone receptor and gene regulated in breast cancer 1 (GREB1) to an equivalent extent as the selective estrogen receptor degrader fulvestrant (Item No. 10011269) in MCF-7 cells when used at concentrations ranging from 1 to 300 nM.

Reference

1. Hu, J., Hu, B., Wang, M., *et al.* Discovery of ERD-308 as a highly potent proteolysis targeting chimera (PROTAC) degrader of estrogen receptor (ER). *J. Med. Chem.* **62**(3), 1420-1442 (2019).

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, 05/09/2024

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