

PRODUCT INFORMATION

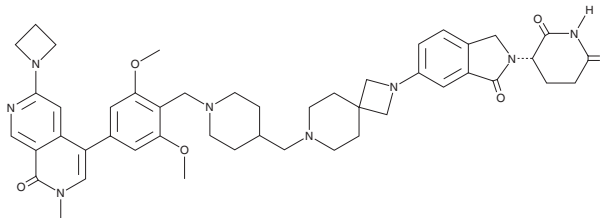


FHD-609

Item No. 46304

CAS Registry No.: 2676211-64-4
Formal Name: (3S)-3-[6-[7-[[1-[[4-[6-(1-azetidiny)-1,2-dihydro-2-methyl-1-oxo-2,7-naphthyridin-4-yl]-2,6-dimethoxyphenyl)methyl]-4-piperidinyl)methyl]-2,7-diazaspiro[3.5]non-2-yl]-1,3-dihydro-1-oxo-2H-isoindol-2-yl]-2,6-piperidinedione

MF: C₄₇H₅₆N₈O₆
FW: 829.0
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

FHD-609 is supplied as a solid. A stock solution may be made by dissolving the FHD-609 in the solvent of choice, which should be purged with an inert gas. FHD-609 is sparingly soluble (1-10 mg/ml) in a 1:1 solution of acetonitrile:water.

Description

FHD-609 is a proteolysis-targeting chimera (PROTAC) and derivative of dBRD9 (Item No. 45720).¹ It degrades bromodomain-containing protein 9 (BRD9) in SYO-1 synovial sarcoma cells.

Reference

1. Ruppel, S.K., Yang, Z., Lowe, J.T., *et al.* Compounds and uses thereof. *Foghorn Therapeutics, Inc.* US2021/0230190A1 (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, 08/13/2026

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