

PRODUCT INFORMATION

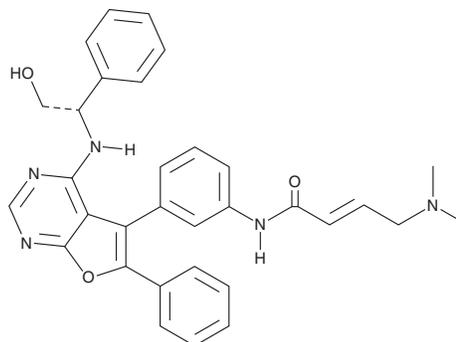


DBPR112

Item No. 44448

CAS Registry No.: 1226549-49-0
Formal Name: (2E)-4-(dimethylamino)-N-[3-[4-[[[(1S)-2-hydroxy-1-phenylethyl]amino]-6-phenylfuro[2,3-d]pyrimidin-5-yl]phenyl]-2-butenamide

Synonym: Gozanertinib
MF: C₃₂H₃₁N₅O₃
FW: 533.6
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

DBPR112 is supplied as a solid. A stock solution may be made by dissolving the DBPR112 in the solvent of choice, which should be purged with an inert gas. DBPR112 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

DBPR112 is an inhibitor of EGFR and dual-mutant EGFR^{L858R/T790M} (IC₅₀s = 15 and 48 nM, respectively).¹ It is selective for these kinases over 384 other kinases but does inhibit 11 additional kinases, including JAK3, Blk, and bone marrow tyrosine kinase on chromosome X (BMX), at 10,000 nM. DBPR112 reduces the viability of HCC827 and H1975 non-small cell lung cancer (NSCLC) cells with 50% cytotoxicity concentration values (CC₅₀s) of 25 and 620 nM, respectively. It decreases tumor weight without affecting body weight in a HCC827 mouse xenograft model when administered at doses of 20 or 50 mg/kg per day.

Reference

1. Lin, S.-Y., Chang Hsu, Y., Peng, Y.-H., *et al.* Discovery of a furanopyrimidine-based epidermal growth factor receptor inhibitor (DBPR112) as a clinical candidate for the treatment of non-small cell lung cancer. *J. Med. Chem.* **62**(22), 10108-10123 (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, 10/28/2025

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