

PRODUCT INFORMATION



JNJ-26483327

Item No. 44468

CAS Registry No.: 807640-87-5
Formal Name: 17-bromo-1,8,9,10,11,12,13,14-octahydro-20-methoxy-13-methyl-4,6-ethenopyrimido[4,5-b][6,1,12]benzoxadiazacyclopentadecine

MF: C₂₂H₂₅BrN₄O₂

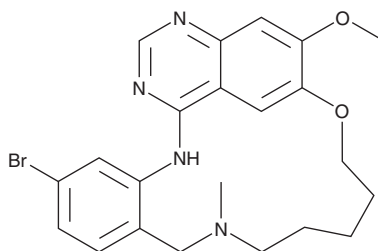
FW: 457.4

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

JNJ-26483327 is supplied as a solid. A stock solution may be made by dissolving the JNJ-26483327 in the solvent of choice, which should be purged with an inert gas. JNJ-26483327 is slightly soluble (0.1-1 mg/ml) in methanol.

Description

JNJ-26483327 is an inhibitor of Met.¹ It also inhibits EGFR, HER2, and HER4 (IC₅₀s = 9.6, 18, and 40.3 nM, respectively). JNJ-26483327 (5 or 10 μM), in combination with trastuzumab, reduces HER2 phosphorylation in BT474 breast cancer cells. It decreases tumor volume when used in combination with trastuzumab in a BT474 mouse xenograft model when administered at a dose of 75 mg/kg twice per day.

Reference

1. Gijzen, M., King, P., Perera, T., *et al.* HER2 phosphorylation is maintained by a PKB negative feedback loop in response to anti-HER2 herceptin in breast cancer. *PLoS Biol.* **8(12)**, e1000563 (2010).

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, 11/04/2025

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