

PRODUCT INFORMATION

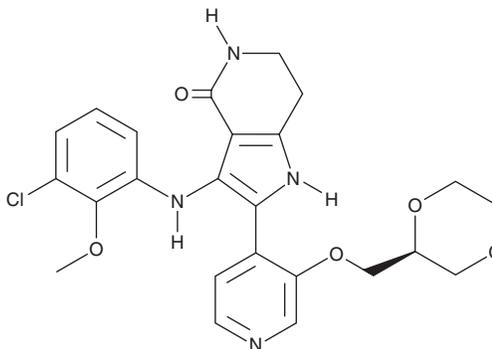


Sevabertinib

Item No. 44586

CAS Registry No.: 2521285-05-0
Formal Name: 3-[[3-chloro-2-methoxyphenyl]amino]-2-[3-[(2S)-1,4-dioxan-2-ylmethoxy]-4-pyridinyl]-1,5,6,7-tetrahydro-4H-pyrrolo[3,2-c]

Synonym: BAY-2927088
MF: C₂₄H₂₅ClN₄O₅
FW: 484.9
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

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

Description

Sevabertinib is a dual inhibitor of wild-type and exon 20 insertion (ex20ins) mutation-containing EGFR and HER2 (IC₅₀s = <0.5 nM for EGFR, EGFR exon20 D770 insSVD, HER2, and HER2 exon20 A775 insYMVA).¹ It is selective for these proteins over HER4 (IC₅₀ = 13.96 nM). Sevabertinib (0.01-1 μM) reduces the viability of A431, NCI-H2170, NCI-H1781, and MDA-MB-175-VII cells endogenously expressing EGFR or HER2 wild-type or insertion mutant gene amplifications. *In vivo*, sevabertinib (5 and 10 mg/kg) reduces tumor growth in a HER2 exon20 A775 insYMVA mutant-containing CTG-2543 patient-derived xenograft (PDX) mouse model. Formulations containing sevabertinib have been used in the treatment of non-small cell lung cancer (NSCLC).

Reference

1. Siegel, F., Siegel, S., Kotýnková, K., *et al.* Sevabertinib, a reversible HER2 inhibitor with activity in lung cancer. *Cancer Discov.* (2025).

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, 12/03/2025

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