

PRODUCT INFORMATION

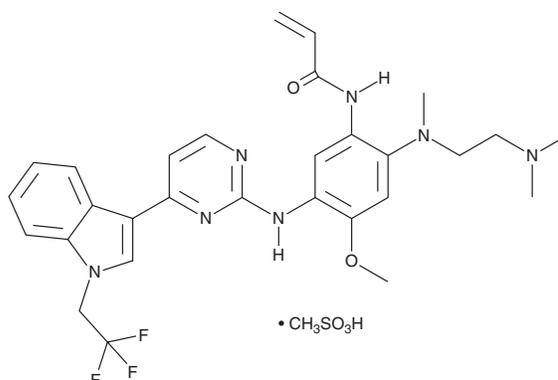


Befotertinib (mesylate)

Item No. 44845

CAS Registry No.: 2226167-02-6
Formal Name: N-[2-[[2-(dimethylamino)ethyl]methylamino]-4-methoxy-5-[[4-[1-(2,2,2-trifluoroethyl)-1H-indol-3-yl]-2-pyrimidinyl]amino]phenyl]-2-propenamide, monomethanesulfonate

Synonym: D-0316
MF: $C_{29}H_{32}F_3N_7O_2 \cdot CH_3SO_3H$
FW: 663.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Befotertinib (mesylate) is supplied as a solid. A stock solution may be made by dissolving the befotertinib (mesylate) in the solvent of choice, which should be purged with an inert gas. Befotertinib (mesylate) is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1). We do not recommend storing the aqueous solution for more than one day.

Description

Befotertinib is an inhibitor of mutant EGFRs.¹ It selectively inhibits EGFR^{L858R}, EGFR^{E709K}, EGFR^{G719A}, and EGFR^{L861Q} (IC_{50} s = 0.2, 14, 1.2, and 13 nM, respectively), as well as EGFR containing an exon 19 deletion and EGFR containing an exon 18 deletion (IC_{50} s = 0.2 and 24 nM, respectively), over wild-type EGFR (IC_{50} = 360 nM).

Reference

1. Oiki, H., Suda, K., Hamada, A., *et al.* Efficacy of conventional and novel tyrosine kinase inhibitors for uncommon EGFR mutations—an in vitro study. *Cells* **14**(17), 1386 (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, 01/26/2026

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