

PRODUCT INFORMATION

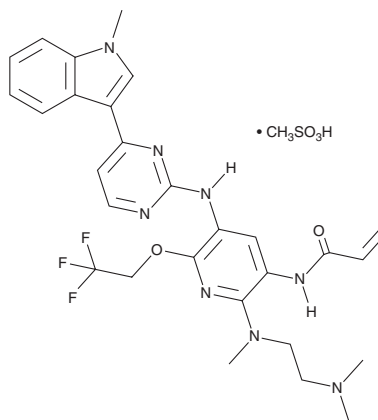


Alflutinib (mesylate)

Item No. 31473

CAS Registry No.: 2130958-55-1
Formal Name: N-[2-[[2-(dimethylamino)ethyl]methylamino]-5-[[4-(1-methyl-1H-indol-3-yl)-2-pyrimidinyl]amino]-6-(2,2,2-trifluoroethoxy)-3-pyridinyl]-2-propenamide, monomethanesulfonate

Synonym: AST2818
MF: C₂₈H₃₁F₃N₈O₂ • CH₃SO₃H
FW: 664.7
Purity: ≥98%
UV/Vis.: λ_{max}: 267, 332 nm
Supplied as: A crystalline 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

Alflutinib (mesylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the alflutinib (mesylate) in the solvent of choice, which should be purged with an inert gas. Alflutinib (mesylate) is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml. Alflutinib (mesylate) is also slightly soluble in DMSO and dimethyl formamide.

Description

Alflutinib is an inhibitor of mutant EGFRs.¹ It is selective for EGFR^{L858R}, EGFR^{G719X}, EGFR^{L858R}, EGFR^{L861A}, and EGFR^{T790M} mutant EGFRs over wild-type EGFR. Alflutinib (10 and 30 mg/kg) reduces tumor growth in an EGFR^{L858R} and EGFR^{T790M}-expressing LU1868 non-small cell lung cancer (NSCLC) patient-derived xenograft (PDX) mouse model.

Reference

1. Shi, Y., Zhang, S., Hu, X., *et al.* Safety, clinical activity, and pharmacokinetics of alflutinib (AST2818) in patients with advanced NSCLC with EGFR T790M mutation. *J. Thorac. Oncol.* **15(6)**, 1015-1026 (2020).

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, 10/18/2022

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