

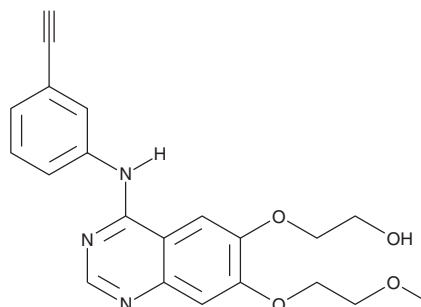
PRODUCT INFORMATION



Desmethyl Erlotinib

Item No. 23394

CAS Registry No.: 183321-86-0
Formal Name: 2-[[4-[(3-ethynylphenyl)amino]-7-(2-methoxyethoxy)-6-quinazolinyloxy]-ethanol
Synonyms: CP 473,420, O-Desmethyl Erlotinib, OSI-420
MF: C₂₁H₂₁N₃O₄
FW: 379.4
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 247, 335 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

Desmethyl erlotinib is supplied as a crystalline solid. A stock solution may be made by dissolving the desmethyl erlotinib in the solvent of choice, which should be purged with an inert gas. Desmethyl erlotinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of desmethyl erlotinib in these solvents is approximately 0.25, 25, and 50 mg/ml, respectively.

Desmethyl erlotinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, desmethyl erlotinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Desmethyl erlotinib has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Desmethyl erlotinib is a metabolite of erlotinib (Item No. 10483).¹

References

1. Sparidans, R.Q., Rosing, H., Rood, J.M., *et al.* Liquid chromatography-tandem mass spectrometric assay for therapeutic drug monitoring of the B-Raf inhibitor encorafenib, the EGFR inhibitors afatinib, erlotinib and gefitinib and the O-desmethyl metabolites of erlotinib and gefitinib in human plasma. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1033-1034**, 390-398 (2016).

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, 03/02/2023

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