

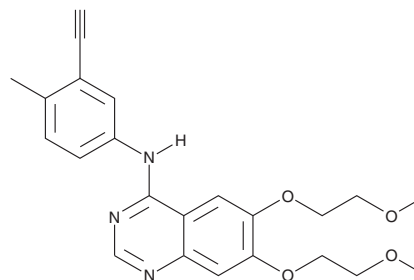
PRODUCT INFORMATION



4-methyl Erlotinib

Item No. 9001510

CAS Registry No.: 1346601-52-2
Formal Name: N-(3-ethynyl-4-methylphenyl)-6,7-bis(2-methoxyethoxy)-4-quinazolinamine
MF: $C_{23}H_{25}N_3O_4$
FW: 407.5
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 226, 248, 336 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

4-methyl Erlotinib is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-methyl erlotinib in the solvent of choice. 4-methyl Erlotinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 4-methyl erlotinib in these solvents is approximately 0.25, 25, and 50 mg/ml, respectively.

4-methyl Erlotinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-methyl erlotinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 4-methyl 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

4-methyl Erlotinib is a methylated derivative of erlotinib (Item Nos. 10483 | 35517).^{1,2} It has been used as an internal standard for the quantification of erlotinib and erlotinib metabolites in human plasma

References

1. Lepper, E.R., Swain, S.M., Tan, A.R., *et al.* Liquid-chromatographic determination of erlotinib (OSI-774), an epidermal growth factor receptor tyrosine kinase inhibitor. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **796**(1), 181-188 (2003).
2. Svedberg, A., Gr  n, H., Vikstr  m, A., *et al.* A validated liquid chromatography tandem mass spectrometry method for quantification of erlotinib, OSI-420 and didesmethyl erlotinib and semi-quantification of erlotinib metabolites in human plasma. *J. Pharm. Biomed. Anal.* **107**, 186-195 (2015).

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, 04/01/2025

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