

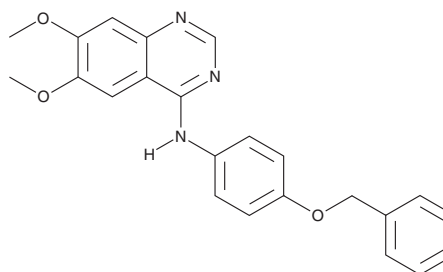
PRODUCT INFORMATION



EGFR/ErbB2 Inhibitor

Item No. 17568

CAS Registry No.: 179248-61-4
Formal Name: 6,7-dimethoxy-N-[4-(phenylmethoxy)phenyl]-4-quinazolinamine
MF: C₂₃H₂₁N₃O₃
FW: 387.4
Purity: ≥95%
UV/Vis.: λ_{max}: 218, 246, 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

EGFR/ErbB2 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the EGFR/ErbB2 inhibitor in the solvent of choice, which should be purged with an inert gas. EGFR/ErbB2 inhibitor is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of EGFR/ErbB2 inhibitor in these solvents is approximately 1 and 10 mg/ml, respectively.

EGFR/ErbB2 inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, EGFR/ErbB2 inhibitor should first be dissolved in DMF and then diluted with the aqueous buffer of choice. EGFR/ErbB2 inhibitor has a solubility of approximately 0.1 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

EGFR/ErbB2 inhibitor is a cell-permeable inhibitor of EGFR and c-ErbB2 (HER2, Neu; IC₅₀s = 20 and 79 nM, respectively).¹ It inhibits the proliferation of cancer cells overexpressing either c-ErbB2 or EGFR (IC₅₀s = 2.3-2.5 μM).¹ EGFR/ErbB2 inhibitor displays a high level of selectivity for these kinases.²

References

1. Cockerill, S., Stubberfield, C., Stables, J., *et al.* Indazolylamino quinazolines and pyridopyrimidines as inhibitors of the EGFR and C-erbB-2. *Bioorg. Med. Chem. Lett.* **11(11)**, 1401-1405 (2001).
2. Anastassiadis, T., Deacon, S.W., Devarajan, K., *et al.* Comprehensive assay of kinase catalytic activity reveals features of kinase inhibitor selectivity. *Nat. Biotechnol.* **29(11)**, 1039-1045 (2011).

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, 01/03/2023

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