

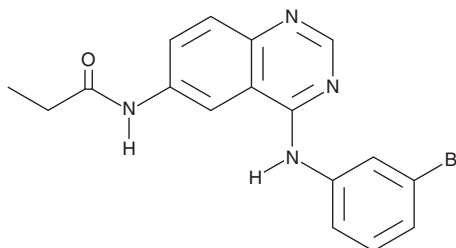
PRODUCT INFORMATION



PD 174265

Item No. 17649

CAS Registry No.: 216163-53-0
Formal Name: N-[4-[(3-bromophenyl)amino]-6-quinazoliny]-propanamide
MF: C₁₇H₁₅BrN₄O
FW: 371.2
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 306, 346 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

PD 174265 is supplied as a crystalline solid. A stock solution may be made by dissolving the PD 174265 in the solvent of choice, which should be purged with an inert gas. PD 174265 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PD 174265 in these solvents is approximately 1, 10, and 30 mg/ml, respectively.

Description

PD 174265 is a potent, cell-permeable inhibitor of the tyrosine kinase activity of the epidermal growth factor (EGF) receptor (IC₅₀ = 0.45 nM).¹ It is effective in cells, blocking tyrosine phosphorylation induced by either EGF or heregulin (IC₅₀s = 39 and 220 nM, respectively).¹ The action of PD 174265 is reversible, which makes it a useful tool for comparative studies with irreversible inhibitors of EGFR kinase activity.¹

Reference

1. Fry, D.W., Bridges, A.J., Denny, W.A., *et al.* Specific, irreversible inactivation of the epidermal growth factor receptor and erbB2, by a new class of tyrosine kinase inhibitor. *Proc. Natl. Acad. Sci. USA* **95**, 12022-12027 (1998).

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, 11/17/2022

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