

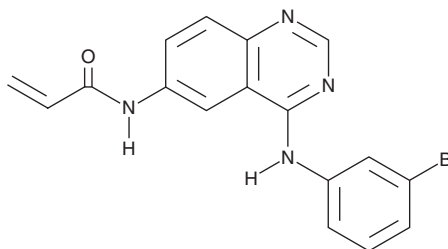
PRODUCT INFORMATION



PD 168393

Item No. 21059

CAS Registry No.: 194423-15-9
Formal Name: N-[4-[(3-bromophenyl)amino]-6-quinazoliny]-2-propenamide
MF: C₁₇H₁₃BrN₄O
FW: 369.2
Purity: ≥95%
UV/Vis.: λ_{max}: 243, 317, 347 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 168393 is supplied as a crystalline solid. A stock solution may be made by dissolving the PD 168393 in the solvent of choice, which should be purged with an inert gas. PD 168393 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PD 168393 in these solvents is approximately 1, 10, and 30 mg/ml, respectively.

Description

PD 168393 is a cell-permeable, irreversible inhibitor of EGFR kinase activity (IC₅₀ = 0.70 nM).¹ It is effective *in vivo*, suppressing the growth of human epidermoid carcinoma xenografts in mice.¹ PD 168393 is used in cells and *in vivo* to irreversibly inhibit EGFR signaling.²⁻⁴

References

1. Fry, D.W., Kraker, A.J., McMichael, A., *et al.* A specific inhibitor of the epidermal growth factor receptor tyrosine kinase. *Science* **265**(5175), 1093-1095 (1994).
2. Pulit-Penaloza, J.A., Sapkota, B., Esser, E.S., *et al.* Modulation of influenza vaccine immune responses using an epidermal growth factor receptor kinase inhibitor. *Sci. Rep.* **5**, 12321 (2015).
3. Sun, X., Liang, J., Yao, X., *et al.* The activation of EGFR promotes myocardial tumor necrosis factor-α production and cardiac failure in endotoxemia. *Oncotarget* **6**(34), 35478-35495 (2015).
4. Wang, S.H., Wang, S.C., Chen, P.C., *et al.* Induction of cyclooxygenase-2 gene by *Candida albicans* through EGFR, ERK, and p38 pathways in human urinary epithelium. *Med. Mycol.* [Epub ahead of print] (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, 11/07/2022

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