

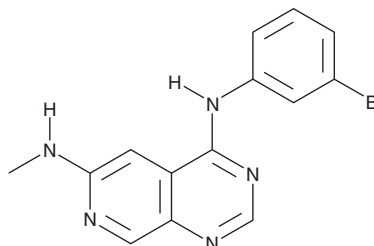
PRODUCT INFORMATION



PD 158780

Item No. 13329

CAS Registry No.: 171179-06-9
Formal Name: N⁴-(3-bromophenyl)-N⁶-methyl-
pyrido[3,4-d]pyrimidine-4,6-diamine
MF: C₁₄H₁₂BrN₅
FW: 330.2
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 312, 404 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 158780 is supplied as a crystalline solid. A stock solution may be made by dissolving the PD 158780 in the solvent of choice, which should be purged with an inert gas. PD 158780 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PD 158780 in these solvents is approximately 2, 10, and 30 mg/ml, respectively.

Description

PD 158780 is a pyridopyrimidine derivative that reversibly inhibits auto and transphosphorylation of all four members of the ErbB receptor tyrosine kinase superfamily: EGFR (IC₅₀ = 8 μM), ErbB2 (IC₅₀ = 49 nM), ErbB3, and ErbB4 (IC₅₀ = 52 nM).¹ It does not inhibit PDGF or FGF-mediated tyrosine phosphorylation.¹ At 0.5 μM, PD 158780 has been shown to induce G₁ cell cycle arrest in MCF10A breast cancer cells.²

References

1. Fry, D.W., Nelson, J.M., Slintak, V., *et al.* Biochemical and antiproliferative properties of 4-[Ar(alk)ylamino] pyridopyrimidines, a new chemical class of potent and specific epidermal growth factor receptor tyrosine kinase inhibitor. *Biochem. Pharmacol.* **54**, 877-887 (1997).
2. Gonzales, A.J. and Fry, D.W. G1 cell cycle arrest due to the inhibition of erbB family receptor tyrosine kinases does not require the retinoblastoma protein. *Exp. Cell Res.* **303**, 56-67 (2005).

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, 12/01/2022

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