

PRODUCT INFORMATION



PD 173952

Item No. 45469

CAS Registry No.: 305820-75-1

Formal Name: 6-(2,6-dichlorophenyl)-8-methyl-2-[[4-(4-morpholinyl)phenyl]amino]pyrido[2,3-d]pyrimidin-7(8H)-one

MF: C₂₄H₂₁Cl₂N₅O₂

FW: 482.4

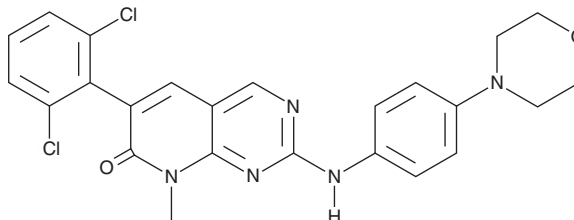
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PD 173952 is supplied as a solid. A stock solution may be made by dissolving the PD 173952 in the solvent of choice, which should be purged with an inert gas. PD 173952 is soluble (≥10 mg/ml) in DMSO.

Description

PD 173952 is a multi-kinase inhibitor.^{1,2} It inhibits LYN, Abl, CSK, and c-Src in cell-free assays (IC₅₀s = 0.3, 1.7, 6.6, and 8.5 nM, respectively). PD 173952 inhibits the growth of HCT-8, SW620, and HT-29 colon carcinoma cells (IC₅₀s = 1.53, 1.8, and 0.526 μM, respectively).¹ It induces cell cycle arrest at the G₂/M phase in HT-29 cells when used at a concentration of 1 μM. PD 173952 (0.5 μM) induces apoptosis in K562 and MEG-01 chronic myeloid leukemia cells.² It also inhibits ristocetin-induced aggregation in human platelet-rich plasma.³

References

1. Kraker, A.J., Hartl, B.G., Amar, A.M., *et al.* Biochemical and cellular effects of c-Src kinase-selective pyrido[2, 3-d]pyrimidine tyrosine kinase inhibitors. *Biochem. Pharmacol.* **60(7)**, 885-898 (2000).
2. Dorsey, J.F., Cunnick, J.M., Lanehart, R., *et al.* Interleukin-3 protects Bcr-Abl-transformed hematopoietic progenitor cells from apoptosis induced by Bcr-Abl tyrosine kinase inhibitors. *Leukemia* **16(9)**, 1589-1595 (2002).
3. Marshall, S.J., Senis, Y.A., Auger, J.M., *et al.* GPIb-dependent platelet activation is dependent on Src kinases but not MAP kinase or cGMP-dependent kinase. *Blood* **103(7)**, 2601-2609 (2004).

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, 03/24/2026

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