PRODUCT INFORMATION



PD 161570

Item No. 24681

CAS Registry No.: Formal Name:	192705-80-9 N-[6-(2,6-dichlorophenyl)-2- [[4-(diethylamino)butyl]amino] pyrido[2,3-d]pyrimidin-7-yl]-N'- (1,1-dimethylethyl)-urea	H I	H ^N O
MF:	C ₂₆ H ₃₅ Cl ₂ N ₇ O	N N	N N H
FW:	532.5	Ĩ	
Purity:	≥98%		Ň
Supplied as:	A solid		
Storage:	-20°C		
Stability:	≥4 years		CI ~ ~

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

Laboratory Procedures

PD 161570 is supplied as a solid. A stock solution may be made by dissolving the PD 161570 in the solvent of choice, which should be purged with an inert gas. PD 161570 is soluble in organic solvents such as DMSO and 0.1 M HCI. The solubility of PD 161570 in these solvents is approximately 100 and 70 mM, respectively.

Description

PD 161570 is an inhibitor of FGF receptor 1 (FGFR1; IC_{50} = 40 nM).¹ It is selective for FGFR1 over PDGF receptor β (PDGF β) and the EGF receptor (EGFR; IC_{50} s = 262 and 3,700 nM, respectively). PD 161570 inhibits constitutive phosphorylation of FGFR1 in Sf9 insect cells overexpressing human FGFR1 and in A121 ovarian carcinoma cells and inhibits growth of A121 cells in a time-dependent manner.

Reference

1. Batley, B.L., Doherty, A.M., Hamby, J.M., et al. Inhibition of FGF-1 receptor tyrosine kinase activity by PD 161570, a new protein-tyrosine kinase inhibitor. Life Sci. 62(2), 143-150 (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.

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, 10/24/2022

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