

PD 173074

Cell-permeable, potent inhibitor of FGF and VEGF receptors.

A cell-permeable compound that acts as a potent, ATP-competitive, and reversible inhibitor of FGF and VEGF receptors. PD 173074 inhibits PDGFR and c-Src only at much higher concentration. Shown to inhibit the autophosphorylation of endogenous FGFR1 and overexpressed VEGFR2.

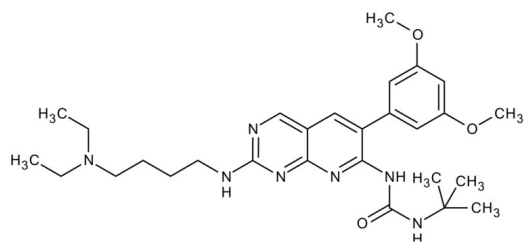
Ordering Information

[Order Online »](#)

ENZ-CHM112-0010	10mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	1-t-Butyl-3-(6-(3,5-dimethoxyphenyl)-2-(4-diethylaminobutylamino)-pyrido[2,3-d]pyrimidin-7-yl)urea
Appearance	White powder solid.
CAS	219580-11-7
Couple Target	FGF receptor, VEGFR
Couple Type	Inhibitor
Formula	$C_{28}H_{41}N_7O_3$
MW	523.7
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com