

PLX4720

Potent and selective B-Raf V600E inhibitor

Equally potent to c-Raf-1 (Y340D and Y341D mutations), and shows 10-fold selectivity for B-Raf V600E over wild-type B-Raf.

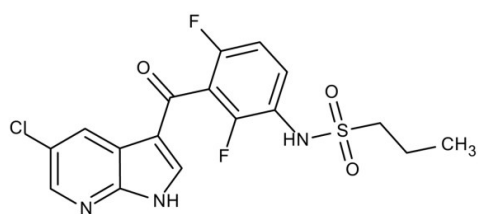
Ordering Information

[Order Online »](#)

ENZ-CHM159-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(3-(5-chloro-1H-pyrrolo[2,3-b]pyridine-3-carbonyl)-2,4-difluorophenyl)propane-1-sulfonamide, N-[3-[(5-Chloro-1H-pyrrolo[2,3-b]pyridin-3-yl)carbonyl]-2,4-difluorophenyl]-1-propanesulfonamide
Appearance	Beige powder.
CAS	918505-84-7
Couple Target	Raf
Couple Type	Inhibitor
Formula	$C_{17}H_{14}ClF_2N_3O_3S$
Identity	Determined by EM-MS, NMR
MW	413.8
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (83mg/ml) or ethanol (1mg/ml).

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-fr@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