

bpV(HOpic)

PTP inhibitor, PTEN Inhibitor

Protein phosphotyrosine phosphatase inhibitor.

Citations: 8

[View Online »](#)

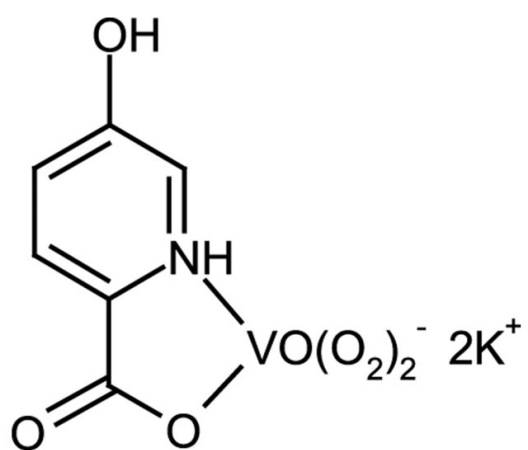
Ordering Information

[Order Online »](#)

ALX-270-206-M005	5mg
ALX-270-206-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow crystalline solid.
Couple Target	PTEN, PTP
Couple Type	Inhibitor
Formula	$K_2[VO(O_2)_2C_6H_4NO_3]$
MW	347.2
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO or water. Note: This product is not stable in water. Starts decomposition immediately after solubilization. Rate of decomposition is slower in dry DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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