

2-Bromoaldisine

Inhibitor of the Raf/MEK-1/MAPK cascade

Inhibitor of the Raf/MEK-1/MAPK cascade (IC_{50} = 539nM).

Citations: 8

[View Online »](#)

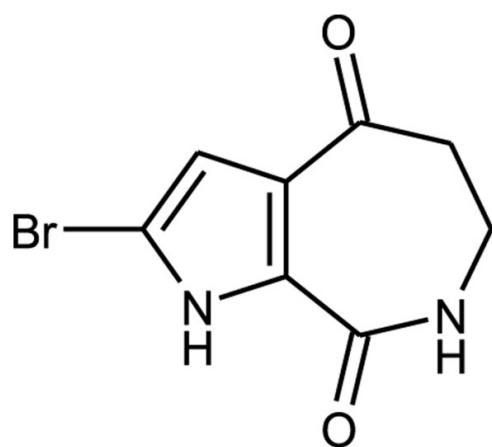
Ordering Information

[Order Online »](#)

ALX-350-388-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	96562-96-8
Couple Target	MAPK
Couple Type	Inhibitor
Formula	$C_8H_7BrN_2O_2$
Identity	Identity determined by 1H-NMR and ESI-MS.
MW	243.1
Purity	≥96% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.
Source	Isolated from sponge <i>Pseudaxynissa</i> sp.

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-
fr@enzolifesciences.com](mailto:info-fr@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)