

Debromohymenialdisine

Chk inhibitor

Inhibitor of G2 DNA damage checkpoint ($IC_{50}=8\mu M$) and check point kinases 1 (Chk1) ($IC_{50}=3\mu M$) and 2 (Chk2) ($IC_{50}=3.5\mu M$). Unlike other checkpoint inhibitors DBH does not inhibit *ataxia-telangiectasia* mutated (ATM) or ATM-Rad3-related protein. Also inhibits MAP kinase kinase 1 (MEK-1) ($IC_{50}=881nM$) but is not as potent as 10Z-hymenialdisine ($IC_{50}=6nM$).

Citations: 3

[View Online »](#)

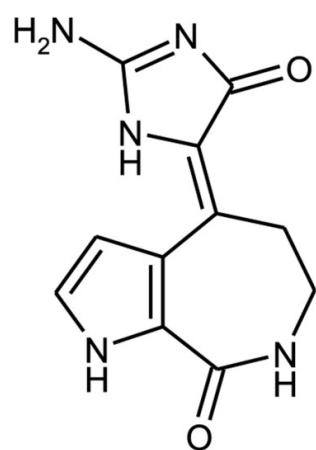
Ordering Information

[Order Online »](#)

ALX-350-290-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DBH, 4-(2-Amino-4-oxo-2-imidazolin-5-ylidene)-4,5,6,7-tetrahydropyrrolo[2,3-c]azepin-8-one
Appearance	Yellow solid.
CAS	75593-17-8
Couple Target	Chk
Couple Type	Inhibitor
Formula	$C_{11}H_{11}N_5O_2$
Identity	Determined by 1H -NMR and MS.
MW	245.2
Purity	≥95% (HPLC)
Solubility	Soluble in 100% ethanol or DMSO.
Source	Isolated from sponge <i>Axinella carteri</i> .

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)