

DMH-1

Selective BMP inhibitor

DMH-1 is an analog of the BMP inhibitor dorsomorphin that potently, selectively inhibits a subset of activin receptor-like kinases, including ALK1, ALK2, ALK3, and ALK6 (IC₅₀ values of 27 nM, 107.9 nM, <5 nM, and 47.6 nM, respectively). It is notably less effective at inhibiting ALK4, ALK5, BMPR2, TGFβR2, AMPK, or KDR (IC50 values of 1825 nM, 565 nM, 3845 nM, 140.4 nM, 1122 nM, and 214.7 nM, respectively). DMH1 is utilized for applications in stem cell biology, specifically for cell differentiation and reprogramming, as it increases cardiomyocyte differentiation in mouse embryonic stem cells (ESCs) and promotes neurogenesis in human-induced pluripotent stem cells (iPSCs).

Ordering Information

[Order Online »](#)

ENZ-CHM324-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 2 years or -20°C for 1 year.
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	BMP Inhibitor II, ALK2/3 Inhibitor, 4-(6-(4-Isopropoxyphenyl)pyrazolo[1,5-a]pyrimidin-3-yl)quinoline
Appearance	Off-white to yellow solid.
CAS	1206711-16-1
Couple Target	Activin receptor
Couple Type	Inhibitor
Formula	$C_{24}H_{20}N_4O$
Identity	Identity determined by LCMS.
MW	380.44
Purity	≥99% (LCMS)
Solubility	Soluble in DMSO (11.5 mg/ml).



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)