# **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 (IC $_{50}$  values of 27 nM, 107.9 nM, <5 nM, and 47.6 nM, respectively). It is notably less effective at inhibiting ALK4, ALK5, BMPR2, TGF $\beta$ R2, 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).

• GMP format available

# **Ordering Information**

**Order Online** »

ENZ-CHM324-0001

1mg

Manuals, SDS & CofA

**View Online »** 

### **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

Inhibitor **Couple Type** 

**Formula**  $C_{24}H_{20}N_4O$ 

Identity Identity determined by LCMS.

380.44 MW

**Purity** ≥99% (LCMS)

**Solubility** Soluble in DMSO (11.5 mg/ml).

eu@enzolifesciences.com