

Dorsomorphin . hydrochloride

Potent and selective **AMPK** inhibitor.

Potent and selective AMPK inhibitor. Dorsomorphin also inhibits the bone morphogenetic protein type 1 receptors ACTR-I (ALK-2), BMPR-IA (ALK-3), and BMPR-IB (ALK-6).

Citations: 4

[View Online »](#)

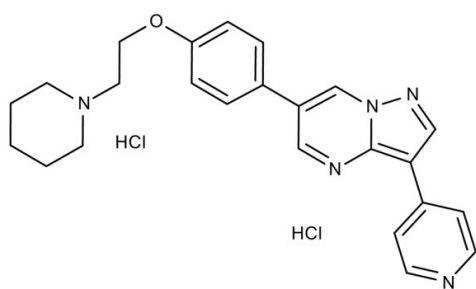
Ordering Information

[Order Online »](#)

ENZ-CHM141-0005	5mg
ENZ-CHM141-0010	10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-[4-[2-(1-Piperidinyloxy)phenyl]-3-(4-pyridinyl)-pyrazolo[1,5-a]pyrimidine dihydrochloride
Appearance	Light yellow powder solid.
CAS	1219168-18-9
Couple Target	Activin receptor, AMPK, BMP receptor
Couple Type	Inhibitor
Formula	$C_{24}H_{25}N_5O \cdot 2 HCl$
MW	472.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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