

DZNep . hydrochloride

EZH2 inhibitor

DZNep hydrochloride is a potent inhibitor of the enzyme EZH2, a histone methyltransferase. This compound acts as an epigenetic modifier by blocking the trimethylation of lysine 27 on histone H3 and lysine 20 on histone H4.

DZNep hydrochloride is also a competitive inhibitor of the Adenosylhomocysteinase (AHCY) S-adenosylhomocysteine hydrolase, with a K_i value of 50 pM. It has shown promising results in inducing selective apoptosis in cancer cells, inhibiting the self-renewal of glioblastoma multiforme cancer stem cells, and enhancing the survival of acute myeloid leukemia blast cells when used in combination with histone deacetylase inhibitors.

Ordering Information

[Order Online »](#)

ENZ-CHM321-0005	5mg
ENZ-CHM321-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. In solvent: Stable at -80°C for 6 months or -20°C for 1 month (sealed and away from moisture).
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial. Sealed storage, avoid moisture.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NSC 617989 . HCl, 3-Deazaneplanocin . HCl
Appearance	White to light brown solid.
CAS	120964-45-6
Couple Target	Adenosylhomocysteinase, Histone methyltransferase
Couple Type	Inhibitor
Formula	$C_{12}H_{15}ClN_4O_3$
MW	298.73
MeltingPoint	168-169°C
Purity	≥99% (HPLC)
Solubility	Soluble in water (50 mg/ml).



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