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 Ki 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 $C_{12}H_{15}CIN_4O_3$

MW 298.73

MeltingPoint 168-169°C

Purity ≥99% (HPLC)

Soluble in water (50 mg/ml).



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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