

3-Deazaneplanocin (DZNep)

Catalog #: 27617-1

Lot #: 230209

Size: 5 mg

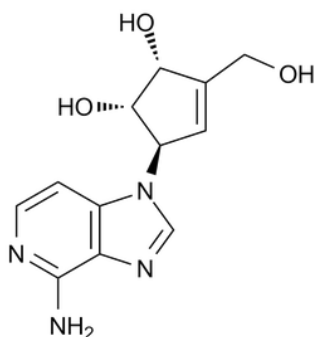
Structure:

CAS Registry #: 102052-95-9

Purity: ≥98%

Chemical Formula: C₁₂H₂₄N₄O₃

Molecular Weight: 262.26 Da



Description: 3-Deazaneplanocin was synthesized as an inhibitor of S-adenosylhomocysteine hydrolase. It is an analog of adenosine and inhibits S-adenosylhomocysteine hydrolase through competing with the substrate, adenosine.

Appearance: A crystalline solid.

Solubility: insoluble in EtOH; ≥17.07 mg/mL in DMSO; ≥17.43 mg/mL in H₂O

Biological Activity: 3-deazaneplanocin A (DZNep), an analog of adenosine, is a competitive inhibitor of S-adenosylhomocysteine hydrolase with K_i of 50 pM.

Storage/Stability: Store at or below -20°C. Stable for at least 1 year when stored as directed. Solutions not stable for long term storage. Avoid freeze/thaw cycles.

Quality Control: The purity was determined by HPLC.

References:

1. Fujiwara T, et al. J Biol Chem. 2014 Mar 21;289(12):8121-34.
2. Cui, B., et al. J Clin Endocrinol Metab. 2014 Jun;99(6):E962-70.
3. Cheng, L.L., et al. Clin Cancer Res. 2012 Aug 1;18(15):4201-12.
4. Nakagawa, S., et al. Oncol Rep. 2014 Feb;31(2):983-8.