

Sinefungin

Catalog No: 14088, 14089

Format: 5 mg, 1 mg

Chemical Properties:

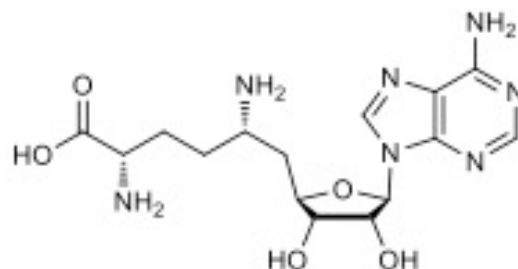
MW = 381.4

C₁₅H₂₃N₇O₅

CAS 58944-73-3

Physical Properties: Tan powder

Names: Sinefungin, 5'-Deoxy-5'-(1,4-diamino-4-carboxybutyl)adenosine; A-9145; Adenosylornithine; Antibiotic 32232RP



Pharmacology: Nucleoside S-adenosyl-1-methionine analog. Potent, competitive methyltransferase (protein, DNA, and RNA methyltransferase) inhibitor (Ref 1,2). IC₅₀ > 1.0 and 2.5 μM for PRMT1 and SET7/9, respectively (ref 3). Binds with greater affinity to the adenine-specific DNA methyltransferase M.TaqI than S-adenosyl-L-homocysteine (ref 4). Inhibits biofilm formation by *Streptococcus pneumoniae* (ref 5).

Solubilization: May be dissolved in water (20 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (HPLC); NMR (Conforms)

References:

1. C Barbes *et al.* *FEMS Microbiol. Lett.* 1990, 57:239
2. MJ Yebra *et al.* *J Antibiot. (Tokyo)* 1991, 44:1141
3. D Cheng *et al.* *J. Biol. Chem.* 2004, 279:23892
4. G Schluckebier *et al.* *J. Mol. Biol.* 1997, 265:56
5. MK Yadav *et al.* *Biomed. Res. Int.* 2014, 2014:156987

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.