

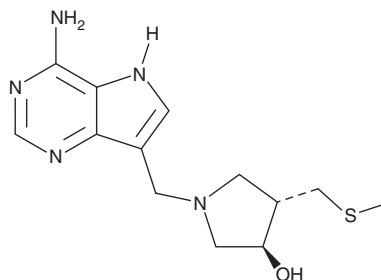
PRODUCT INFORMATION



MT-DADMe-ImmA

Item No. 29711

CAS Registry No.: 653592-04-2
Formal Name: (3R,4S)-1-[(4-amino-5H-pyrrolo[3,2-d]pyrimidin-7-yl)methyl]-4-[(methylthio)methyl]-3-pyrrolidinol
Synonyms: Methylthio-DADMe-Immucillin A, MTDIA
MF: C₁₃H₁₉N₅OS
FW: 293.4
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 279 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MT-DADMe-ImmA is supplied as a crystalline solid. A stock solution may be made by dissolving the MT-DADMe-ImmA in the solvent of choice, which should be purged with an inert gas. MT-DADMe-ImmA is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MT-DADMe-ImmA in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MT-DADMe-ImmA can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of MT-DADMe-ImmA in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MT-DADMe-ImmA is an inhibitor of 5'-methylthioadenosine phosphorylase (MTAP; K_i = 1.7 nM).¹ It increases levels of MTA (Item No. 15593), a product of polyamine metabolism, in FaDu human squamous cell carcinoma cells when used at a concentration of 1 μM.² MT-DADMe-ImmA (5, 9, and 21 mg/kg per day for 28 days) reduces tumor growth in a FaDu mouse xenograft model.

References

1. Evans, G.B., Furneaux, R.H., Lenz, D.H., *et al.* Second generation transition state analogue inhibitors of human 5'-methylthioadenosine phosphorylase. *J. Med. Chem.* **48**(14), 4679-4689 (2005).
2. Basu, I., Cordovano, G., Das, I., *et al.* A transition state analogue of 5'-methylthioadenosine phosphorylase induces apoptosis in head and neck cancers. *J. Biol. Chem.* **282**(29), 21477-21486 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM