

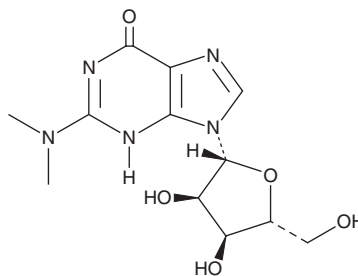
PRODUCT INFORMATION



N²,N²-Dimethylguanosine

Item No. 35347

CAS Registry No.: 2140-67-2
Synonym: N,N-Dimethylguanosine
MF: C₁₂H₁₇N₅O₅
FW: 311.3
Purity: ≥95%
UV/Vis.: λ_{max}: 260 nm
Supplied as: A 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

N²,N²-Dimethylguanosine is supplied as a solid. A stock solution may be made by dissolving the N²,N²-dimethylguanosine in the solvent of choice, which should be purged with an inert gas. N²,N²-Dimethylguanosine is slightly soluble in DMSO.

N²,N²-Dimethylguanosine is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

N²,N²-Dimethylguanosine is a methylated purine nucleoside formed during the degradation of tRNA.¹ Serum levels of N²,N²-dimethylguanosine are elevated in patients with acute myelomonocytic leukemia (AMML) or large cell lung carcinoma.

Reference

1. Mitchell, E.P., Evans, L., Schultz, P., *et al.* Modified nucleosides in human serum. *J. Chromatogr.* **581**(1), 31-40 (1992).

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, 12/12/2022

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