

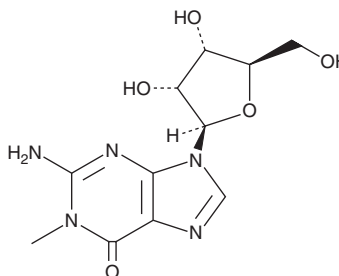
PRODUCT INFORMATION



N¹-Methylguanosine

Item No. 31737

CAS Registry No.: 2140-65-0
Formal Name: 1-methyl-guanosine
Synonyms: 1-Methylguanosine, m¹G, NSC 70897
MF: C₁₁H₁₅N₅O₅
FW: 297.3
Purity: ≥95%
UV/Vis.: λ_{max}: 257 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

N¹-Methylguanosine (m¹G) is supplied as a crystalline solid. A stock solution may be made by dissolving the m¹G in the solvent of choice, which should be purged with an inert gas. m¹G is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of m¹G in these solvents is approximately 1 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 m¹G can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of m¹G in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

m¹G is a methylated purine nucleoside formed during the degradation of tRNA and a biological end product.^{1,2} Levels of m¹G are increased in the urine of patients with malignant tumors compared to those with benign or no tumors and have been used as biomarkers of cancer.^{1,3}

References

1. Liebig, H.M., Müller-Hagedorn, S., Klaus, F., *et al.* Chromatographic, capillary electrophoretic and matrix-assisted laser desorption ionization time-of-flight mass spectrometry analysis of urinary modified nucleosides as tumor markers. *J. Chromatogr. A.* **1071(1-2)**, 271-275 (2005).
2. Mitchell, E.P., Evans, L., Schultz, P., *et al.* Modified nucleosides in human serum. *J. Chromatogr.* **581(1)**, 31-40 (1992).
3. Seidel, A., Brunner, S., Seidel, P., *et al.* Modified nucleosides: An accurate tumour marker for clinical diagnosis of cancer, early detection and therapy control. *Br. J. Cancer.* **94(11)**, 1726-1733 (2006).

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, 12/13/2022

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