

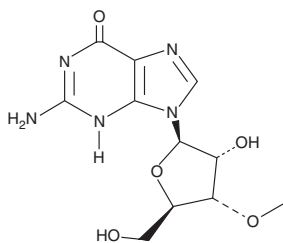
PRODUCT INFORMATION



3'-O-Methylguanosine

Item No. 40292

CAS Registry No.: 10300-27-3
Synonym: NSC 106541
MF: $C_{11}H_{15}N_5O_5$
FW: 297.3
Purity: $\geq 98\%$
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

3'-O-Methylguanosine is supplied as a solid. A stock solution may be made by dissolving the 3'-O-methylguanosine in the solvent of choice, which should be purged with an inert gas. 3'-O-Methylguanosine is soluble in methanol.

Description

3'-O-Methylguanosine is an antiviral nucleoside analog.¹ It inhibits viral plaque formation and viral RNA synthesis in vaccinia virus-infected Vero cells when used at a concentration of 10 μM .

Reference

1. Goswami, B.B. and Sharma, O.K. Inhibition of vaccinia virus growth and virus-specific RNA synthesis by 3'-O-methyl adenosine and 3'-O-methyl guanosine. *J. Virol.* **45**(3), 1164-1167 (1983).

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, 03/19/2024

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