

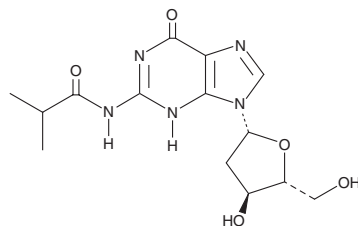
PRODUCT INFORMATION



N²-Isobutyryl-2'-deoxyguanosine

Item No. 40284

CAS Registry No.: 68892-42-2
Formal Name: 2'-deoxy-N-(2-methyl-1-oxopropyl)-guanosine
MF: C₁₄H₁₉N₅O₅
FW: 337.3
Purity: ≥95%
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-Isobutyryl-2'-deoxyguanosine is supplied as a solid. A stock solution may be made by dissolving the N²-isobutyryl-2'-deoxyguanosine in the solvent of choice, which should be purged with an inert gas. N²-Isobutyryl-2'-deoxyguanosine is sparingly soluble (1-10 mg/ml) in ethanol.

Description

N²-Isobutyryl-2'-deoxyguanosine is a nucleoside derivative and synthetic intermediate that has been used in the synthesis of the ppGpp derivative relacin, an inhibitor of *E. coli* RelA with antibacterial activity, as well as nucleoside derivatives with antiviral activity against HIV-1.^{1,2}

References

1. Wexselblatt, E., Oppenheimer-Shaanan, Y., Kaspary, I., *et al.* Relacin, a novel antibacterial agent targeting the Stringent Response. *PLoS Pathog.* **8**(9), e1002925 (2012).
2. Zhang, H.-W., Coats, S.J., Bondada, L., *et al.* Synthesis and evaluation of 3'-azido-2',3'-dideoxypurine nucleosides as inhibitors of human immunodeficiency virus. *Bioorg. Med. Chem. Lett.* **20**, 60-64 (2010).

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, 08/07/2024

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