

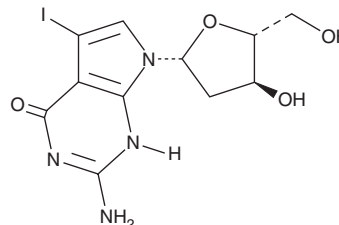
PRODUCT INFORMATION



7-Deaza-7-iodo-2'-deoxyguanosine

Item No. 40643

CAS Registry No.: 172163-62-1
Formal Name: 2-amino-7-(2-deoxy-β-D-erythro-pentofuranosyl)-1,7-dihydro-5-iodo-4H-pyrrolo[2,3-d]pyrimidin-4-one
Synonym: 7-Deaza-2'-deoxy-7-iodoguanosine
MF: C₁₁H₁₃IN₄O₄
FW: 392.2
Purity: ≥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

7-Deaza-7-iodo-2'-deoxyguanosine is supplied as a solid. A stock solution may be made by dissolving the 7-deaza-7-iodo-2'-deoxyguanosine in the solvent of choice, which should be purged with an inert gas. 7-Deaza-7-iodo-2'-deoxyguanosine is slightly soluble (0.1-1 mg/ml) in DMSO

Description

7-Deaza-7-iodo-2'-deoxyguanosine is a derivative of the purine nucleoside 2'-deoxyguanosine (Item No. 9002864). It has been used as an intermediate in the synthesis of fluorescent nucleotide derivatives for incorporation into single- or double-stranded oligonucleotides.¹

Reference

1. Ingale, S.A. and Seela, F. Nucleoside and oligonucleotide pyrene conjugates with 1,2,3-triazolyl or ethynyl linkers: Synthesis, duplex stability, and fluorescence changes generated by the DNA-dye connector. *Tetrahedron* **70(2)**, 380-391 (2014).

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, 03/28/2024

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