

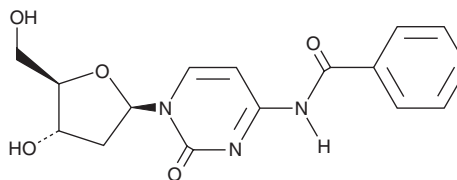
# PRODUCT INFORMATION



## N<sup>4</sup>-benzoyl-2'-Deoxycytidine

Item No. 39934

CAS Registry No.: 4836-13-9  
Formal Name: N-benzoyl-2'-deoxy-cytidine  
Synonym: N<sup>4</sup>-benzoyl-dC  
MF: C<sub>16</sub>H<sub>17</sub>N<sub>3</sub>O<sub>5</sub>  
FW: 331.3  
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

N<sup>4</sup>-benzoyl-2'-Deoxycytidine is supplied as a solid. A stock solution may be made by dissolving the N<sup>4</sup>-benzoyl-2'-deoxycytidine in the solvent of choice, which should be purged with an inert gas. N<sup>4</sup>-benzoyl-2'-Deoxycytidine is soluble in DMSO and methanol.

### Description

N<sup>4</sup>-benzoyl-2'-Deoxycytidine is a building block and derivative of the pyrimidine nucleoside 2'-deoxycytidine (Item Nos. 34708 | 30125).<sup>1,2</sup> It has been used in the synthesis of nucleosides and nucleotides with anticancer activity, as well as oligonucleotides.

### References

1. Yamaguchi, T. and Saneyoshi, M. Synthetic nucleosides and nucleotides. XXI. On the synthesis and biological evaluations of 2'-deoxy- $\alpha$ -D-ribofuranosyl nucleosides and nucleotides. *Chem. Pharm. Bull. (Tokyo)* **32(4)**, 1441-1450 (1984).
2. Cazenave, C., Bathany, K., and Rayner, B. Formation of N-branched oligonucleotides as by-products in solid-phase oligonucleotide synthesis. *Oligonucleotides* **16(2)**, 181-185 (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, 03/08/2024

#### 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