

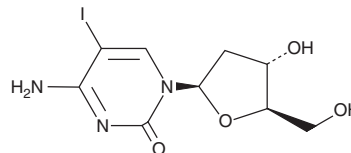
# PRODUCT INFORMATION



## 5-Iodo-2'-deoxycytidine

Item No. 35696

**CAS Registry No.:** 611-53-0  
**Formal Name:** 2'-deoxy-5-iodo-cytidine  
**Synonyms:** ICDR, IdCyd, 5-Iododeoxycytidine, NSC 527083  
**MF:** C<sub>9</sub>H<sub>12</sub>I N<sub>3</sub>O<sub>4</sub>  
**FW:** 353.1  
**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

5-Iodo-2'-deoxycytidine is supplied as a solid. A stock solution may be made by dissolving the 5-iodo-2'-deoxycytidine in the solvent of choice, which should be purged with an inert gas. 5-Iodo-2'-deoxycytidine is slightly soluble in acetonitrile.

5-Iodo-2'-deoxycytidine is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

5-Iodo-2'-deoxycytidine is a nucleoside analog.<sup>1</sup> It inhibits the growth of P815 mastocytoma cells (IC<sub>50</sub> = 14 μM), as well as sensitizes P815 cells to irradiation when used at a concentration of 1 μM.

### Reference

1. Cramer, J.W., Prusoff, W.H., Welch, A.D., *et al.* Studies on the biochemical pharmacology of 5-iodo-2'-deoxycytidine *in vitro* and *in vivo*. *Biochem. Pharmacol.* **11**, 761-768 (1962).

#### 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, 08/20/2024

#### 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