

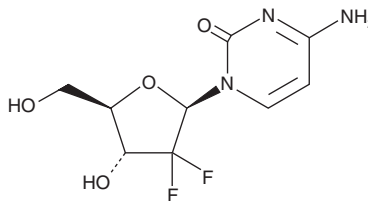
# PRODUCT INFORMATION



## Gemcitabine

Item No. 11690

CAS Registry No.:	95058-81-4
Formal Name:	2'-deoxy-2',2'-difluoro-cytidine
Synonyms:	DDFC, LY 188011, NSC 613327
MF:	C <sub>9</sub> H <sub>11</sub> N <sub>3</sub> O <sub>4</sub> F <sub>2</sub>
FW:	263.2
Purity:	≥98%
UV/Vis.:	λ <sub>max</sub> : 243, 269 nm
Supplied as:	A crystalline solid
Storage:	-20°C
Stability:	≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Gemcitabine is supplied as a crystalline solid. A stock solution may be made by dissolving the gemcitabine in the solvent of choice. Gemcitabine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of gemcitabine in these solvents is approximately 0.25, 5, and 0.1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of gemcitabine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of gemcitabine in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Gemcitabine is an anticancer nucleoside analog that inhibits the growth of HL-60 promyelocytic leukemia cells with an LC<sub>50</sub> value of 40 nM.<sup>1</sup> It inhibits the growth of MX-1 mammary, CX-1, HC-1, GC3, and VRC5 colon, LX-1, Calu-6, and NCI-H460 lung, and HS766T, PaCa-2, PANC-1, and BxPC-3 pancreatic cancer tumors in mouse xenograft models (45-93% inhibition).<sup>2</sup> Gemcitabine is a prodrug that is metabolized to a diphosphate and triphosphate form in cells. The triphosphate form is incorporated into DNA which induces masked chain termination and cell death.<sup>3</sup> By specifically inhibiting growth arrest and DNA damage inducible protein 45 a (Gadd45a), a key mediator of active DNA demethylation, gemcitabine, at concentrations ranging from 34 to 134 nM, inhibits repair-mediated DNA demethylation in a methylation-sensitive reporter assay. Gemcitabine also has broad antiretroviral activity, decreasing MuLV cell infectivity, a murine AIDS model, in cell culture (EC<sub>50</sub> = ~1.5 nM) and inhibits the progression of murine AIDS *in vivo* at a dose of 1-2 mg/kg per day.<sup>4</sup>

### References

1. Ross, D.D. and Cuddy, D.P. *Biochem. Pharmacol.* **48(8)**, 1619-1630 (1994).
2. Merriman, R.L., Hertel, L.W., Schultz, R.M., *et al. Invest. New. Drugs* **14(3)**, 243-247 (1996).
3. Schäfer, A., Schomacher, L., Barreto, G., *et al. PLoS One* **5(11)**, 1-9 (2010).
4. Clouser, C.L., Holtz, C.M., Mullett, M., *et al. PLoS One* **6(1)**, 1-8 (2011).

#### 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, 12/21/2017

#### 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