

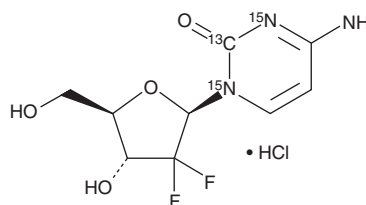
# PRODUCT INFORMATION



## Gemcitabine-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> (hydrochloride)

Item No. 22560

**CAS Registry No.:** 2757566-59-7  
**Formal Name:** 2'-deoxy-2',2'-difluoro-cytidine-2-<sup>13</sup>C-1,3-<sup>15</sup>N<sub>2</sub>, monohydrochloride  
**Synonym:** dFdC  
**MF:** C<sub>8</sub>[<sup>13</sup>C]H<sub>11</sub>F<sub>2</sub>N[<sup>15</sup>N<sub>2</sub>]O<sub>4</sub> • HCl  
**FW:** 302.6  
**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

Gemcitabine-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the gemcitabine-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> (hydrochloride) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Gemcitabine-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> is intended for use as an internal standard for the quantification of gemcitabine (Item Nos. 11690 | 9003096) by GC- or LC-MS. Gemcitabine is an active metabolite of the gemcitabine prodrugs NUC-1031 (Item No. 9003247) and gemcitabine elaidate (Item No. 28303) and a prodrug form of gemcitabine di- and triphosphates.<sup>1-3</sup> Gemcitabine is phosphorylated by intracellular kinases to the intermediate metabolite gemcitabine monophosphate (Item No. 31726) and the active di- and triphosphate forms.<sup>3</sup> It is cytotoxic to HepG2 hepatocellular carcinoma and A549 non-small cell lung cancer (NSCLC) cells (IC<sub>50</sub>s = 5.2 and 16 nM, respectively) and inhibits tumor growth in various breast, colon, lung, and pancreatic cancer mouse xenograft models.<sup>4</sup> Gemcitabine (12 mg/kg) sensitizes tumors to antibodies targeting programmed cell death protein 1 (PD-1), decreases the number of tumor-infiltrating regulatory T cells (Tregs), and increases survival in an MC-38 syngeneic mouse model of colon carcinoma.<sup>5</sup> Formulations containing gemcitabine have been used in the treatment of cancer.

### References

1. Slusarczyk, M., Lopez, M.H., Balzarini, J., *et al.* Application of ProTide technology to gemcitabine: A successful approach to overcome the key cancer resistance mechanisms leads to a new agent (NUC-1031) in clinical development. *J. Med. Chem.* **57**(4), 1531-1542 (2014).
2. Bergman, A.M., Adema, A.D., Balzarini, J., *et al.* Antiproliferative activity, mechanism of action and oral antitumor activity of CP-4126, a fatty acid derivative of gemcitabine, in in vitro and in vivo tumor models. *Invest. New Drugs* **29**(3), 456-466 (2011).
3. Veltkamp, S.A., Pluim, D., van Eijndhoven, M.A.J., *et al.* New insights into the pharmacology and cytotoxicity of gemcitabine and 2',2'-difluorodeoxyuridine. *Mol. Cancer Ther.* **7**(8), 2415-2425 (2008).
4. Merriman, R.L., Hertel, L.W., Schultz, R.M., *et al.* Comparison of the antitumor activity of gemcitabine and ara-C in a panel of human breast, colon, lung and pancreatic xenograft models. *Invest. New. Drugs* **14**(3), 243-247 (1996).
5. Obradovic, A., Ager, C., Turunen, M., *et al.* Systematic elucidation and pharmacological targeting of tumor-infiltrating regulatory T cell master regulators. *Cancer Cell* **41**(5), 933-949 (2023).

#### 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, 05/17/2023

#### 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