

Gemcitabine . hydrochloride

DNA synthesis inhibitor

Deoxycytidine analog with cytotoxic effects. Induces apoptosis and inhibits DNA synthesis.

Citations: 6

[View Online »](#)

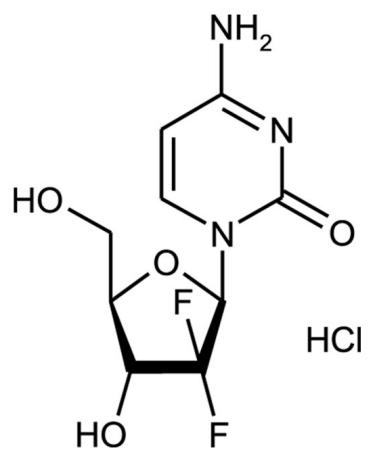
Ordering Information

[Order Online »](#)

ALX-480-101-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2'-Deoxy-2',2'-difluorocytidine . HCl, LY-188011 . HCl
Appearance	White to off-white crystalline powder.
CAS	122111-03-9
Formula	$C_9H_{11}F_2N_3O_4 \cdot HCl$
Identity	Determined by IR.
MI	14: 4386
MW	263.2 . 36.5
Purity	≥97.5% (Assay)
RTECS	HA3840000
Solubility	Soluble in water; slightly soluble in methanol; almost insoluble in 100% ethanol or polar organic solvents.
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)