

Fludarabine

F-ara-A precursor

Prodrug converted to the free nucleoside 9- β -D-arabinosyl-2-fluoroadenine (F-ara-A). F-ara-A enters cells and accumulates as the 5'-triphosphate. Interferes with DNA synthesis and repair. Induces apoptosis of cancer cells. Inhibits RNA transcription.

Citations: 9

[View Online »](#)

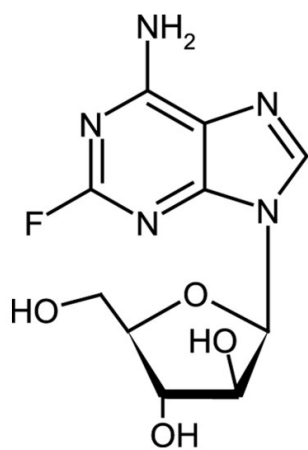
Ordering Information

[Order Online »](#)

ALX-480-100-M005	5mg
ALX-480-100-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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-Fluoroadenine-9-β-D-arabinofuranoside

Appearance White to pale yellow crystalline powder.

CAS 21679-14-1

Formula $C_{10}H_{12}FN_5O_4$

Identity Determined by IR.

MI **14:** 4126

MW 285.2

Purity ≥98% (HPLC)

RTECS AU6207000

Solubility Sparingly soluble in water or organic solvents.

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-
fr@enzolifesciences.com](mailto:info-fr@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)