

DL- α - Difluoromethylornithine . hydrochloride . monohydrate

Inhibitor of ornithine decarboxylase and arginase
Specific irreversible inhibitor of ornithine decarboxylase (ODC), the rate-limiting enzyme in polyamine biosynthesis. Potent chemopreventive agent. Induces apoptosis. Arginase activity inhibitor.

Citations: 15

[View Online »](#)

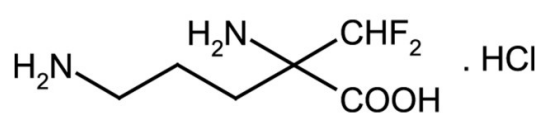
Ordering Information

[Order Online »](#)

ALX-270-283-M010	10mg
ALX-270-283-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DFMO . HCl . H ₂ O, Eflornithine . HCl . H ₂ O, RMI-71782
Appearance	White to off-white solid.
CAS	96020-91-6
Couple Target	Arginase, Ornithine decarboxylase
Couple Type	Inhibitor
Formula	$C_6H_{12}F_2N_2O_2 \cdot HCl \cdot H_2O$
MI	14: 3522
MW	236.6
Purity	≥98% (TLC)
Solubility	Soluble in water (25mg/ml).

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)