

Nutlin-3

Inhibitor of p53/MDM2 interaction

(+/-) -Nutlin-3 is an inhibitor of p53/MDM2 interaction ($IC_{50}=0.09\mu M$).

Citations: 8

[View Online »](#)

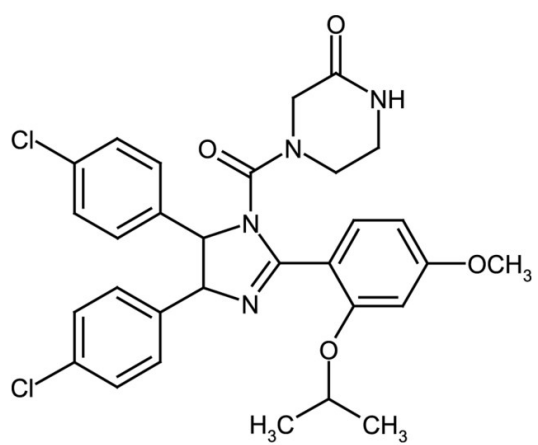
Ordering Information

[Order Online »](#)

ALX-430-128-M001	1mg
ALX-430-128-M005	5mg
ALX-430-128-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored at -20°C. We do not recommend storing the aqueous solution for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-4-[4,5-bis-(4-Chlorophenyl)-2-(2-isopropoxy-4-methoxyphenyl)-4,5-dihydroimidazole-1-carbonyl]-piperazin-2-one
Appearance	White to off-white crystalline solid.
CAS	548472-68-0
Formula	$C_{30}H_{30}Cl_2N_4O_4$
MW	581.5
Purity	≥98% (HPLC (UV))
Solubility	Soluble in methanol, 100% ethanol, DMSO, dimethyl formamide or acetonitrile; sparingly soluble in a 1:10 solution of ethanol:PBS, pH 7.2.

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)