

AM 281

CB₁ receptor ligand

Analog of the cannabinoid receptor (CB) antagonist SR 141716A. Potent CB₁ receptor antagonist/inverse agonist ($K_i=14\text{nM}$).

Citations: 6

[View Online »](#)

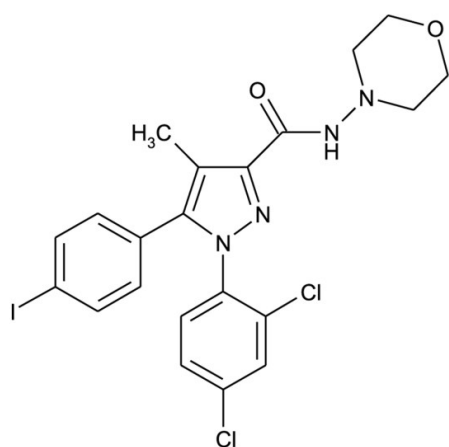
Ordering Information

[Order Online »](#)

ALX-270-240-M001	1mg
ALX-270-240-M005	5mg

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 1-(2,4-Dichlorophenyl)-5-(4-iodophenyl)-4-methyl-N-4-morpholinyl-1H-pyrazole-3-carboxamide

Appearance White to off-white solid.

CAS 202463-68-1

Couple Target Cannabinoid receptor

Couple Type Ligand

Formula $C_{21}H_{19}Cl_2IN_4O_2$

Identity Identity determined by 1H-NMR.

MW 557.2

Purity ≥98% (HPLC)

Solubility Soluble in methanol (1mg/ml) or DMSO (2mg/ml at RT, 50mg/ml in hot DMSO, after cooling this solution is stable over several days.)

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)