

AM-630

Cannabinoid receptor ligand

Acts as a cannabinoid receptor antagonist in mouse brain, vas deferens and guinea pig brain, but an agonist in guinea pig ileum. Acts as an inverse agonist on cloned human CB₁ receptors.

Citations: 7

[View Online »](#)

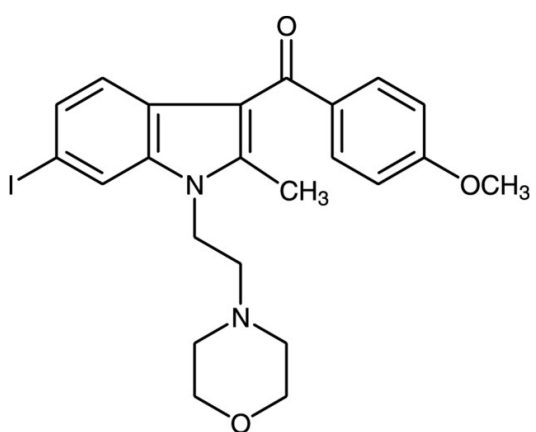
Ordering Information

[Order Online »](#)

BML-CR107-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Store under inert gas and avoid exposure to light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

CAS 164178-33-0

Couple Target Cannabinoid receptor

Couple Type Ligand

Formula $C_{23}H_{25}IN_2O_3$

MW 504.4

Purity ≥98% (TLC)

Solubility Soluble in DMSO or dimethyl formamide.

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)