

AM-251

Cannabinoid receptor ligand

Cannabinoid CB₁ receptor antagonist ($K_i=7.5$ nM). Displays a high level of selectivity (306-fold) for CB₁ over CB₂ receptors. Structurally related to the cannabinoid receptor (CB) antagonist SR 141716A.

Citations: 10

[View Online »](#)

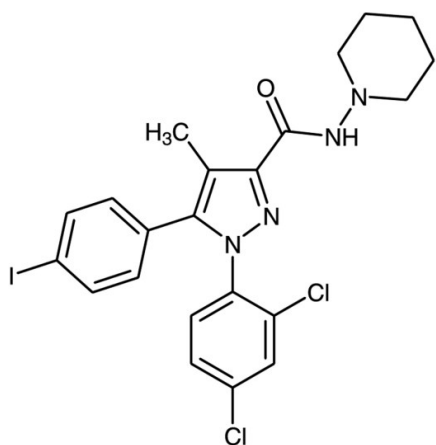
Ordering Information

[Order Online »](#)

BML-CR108-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to Off White
CAS	183232-66-8
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{22}H_{21}Cl_2IN_4O$
Identity	Determined by 1H -NMR.
MW	555.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol 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)