

CP-55,940

Cannabinoid receptor ligand

Cannabinoid CB₁ and CB₂ agonist which is more potent than δ 9-THC. K_d's for cloned/transfected CB₁ and CB₂: 2.6 & 3.7 nM. Has been reported to induce the immediate early gene Krox-24 in corpus striatum and cultured astrocytes. Stimulates proliferation of stimulated B-cells at low nM concentrations.

Citations: 7

[View Online »](#)

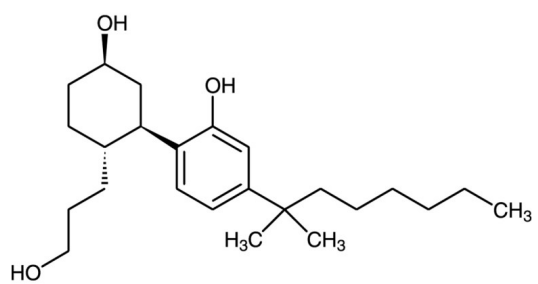
Ordering Information

[Order Online »](#)

BML-CR102-0005	5mg
BML-CR102-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	83002-04-4
Couple Target	Cannabinoid receptor
Couple Type	Ligand
Formula	$C_{24}H_{40}O_3$
MW	376.6
Purity	≥96% (HPLC)
Solubility	Soluble in DMSO (4mg/ml) or ethanol.



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-
be@enzolifesciences.com](mailto:info-be@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)