

# WIN-55,212-2

## Cannabinoid receptor agonist

High affinity cannabinoid receptor (CB) agonist ( $K_i=62.3\text{nM}$  and  $3.3\text{nM}$  at the human cloned  $\text{CB}_1$  and  $\text{CB}_2$  receptors, respectively). Potent analgesic in a rat model of neuropathic pain. Activates p42 and p44 MAP kinase *via* a receptor-mediated signaling.

Citations: 13

[View Online »](#)

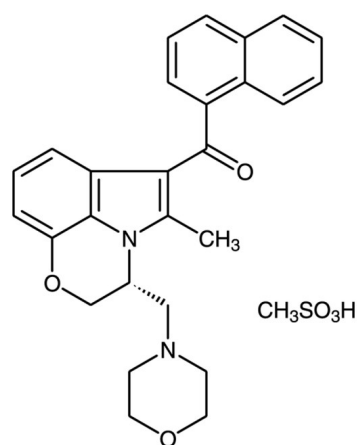
## Ordering Information

[Order Online »](#)

BML-CR105-0010	10mg
BML-CR105-0050	50mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	(R)-(+)-[2,3-Dihydro-5-methyl-3-(4-morpholinylmethyl)pyrrolo[1,2,3-de]-1,4-benzoxazin-6-yl]-1-naphthalenylmethanone . mesylate
Appearance	White to off-white solid.
CAS	131543-23-2
Couple Target	Cannabinoid receptor, MAPK
Couple Type	Activator, Ligand
Formula	$C_{27}H_{26}N_2O_3 \cdot CH_3SO_3H$
MW	522.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO; insoluble in water.

## Technical Info / Product Notes

WIN-55,212-2 is a controlled substance in some countries and may require an import/export license from the local authorities depending on the country of import. Orders for this product will incur an administrative fee to cover license expenses. Please inquire for additional information.



ENZO LIFE SCIENCES,  
INC.  
May 29, 2024  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)