WIN-55,212-2

Cannabinoid receptor agonist

High affinity cannabinoid receptor (CB) agonist (K_i =62.3nM and 3.3nM at the human cloned CB₁ and CB₂ receptors, respectively). Potent analgesic in a rat model of neuropathic pain. Activates p42 and p44 MAP kinase *via* a receptor-mediated signaling.

Citations: 12

View Online »

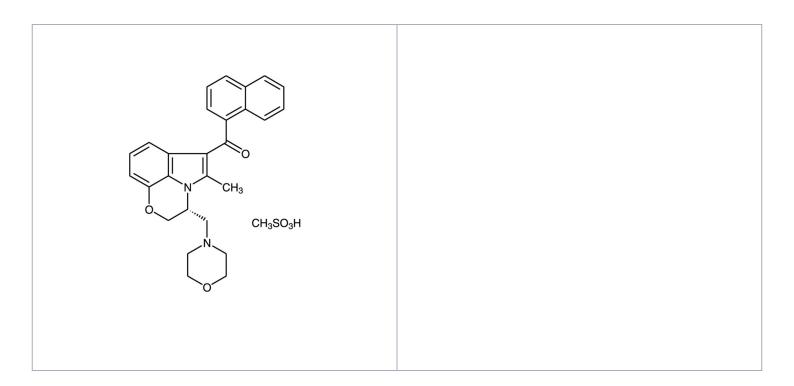
Ordering Information

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 $C_{27}H_{26}N_2O_3$. CH_3SO_3H

MW 522.6

Purity ≥98% (HPLC)

Solubility Soluble in DMSO; insoluble in water.

Technical Info / Product NotesWIN-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.

