

Salvinorin A

κ-opioid receptor ligand

Potent naturally occurring diterpenoid κ1-opioid selective agonist. Displays high affinity at κ1-opioid receptors ($K_i=4.3$ nM).

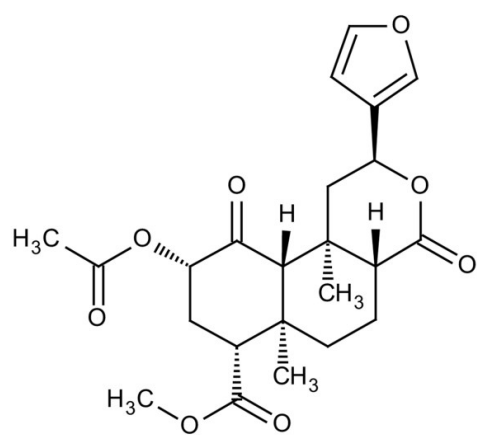
Ordering Information

[Order Online »](#)

BML-NO109-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stored at -20°C, protected from light and oxygen: dry powder is stable for years, solutions are stable up to 2 days (light protected).
Short Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White powder.
CAS	83729-01-5
Couple Target	Opioid receptor
Couple Type	Ligand
MW	432.5
Purity	≥98% (HPLC)
Solubility	Soluble in methanol, pyridine, acetonitrile, and chloroform.
Source	Isolated from cultivated plant <i>Salvia divinorum</i> .

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-
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)