

Dihydroergocristine . mesylate

Serotonin receptor ligand

Serotonin antagonist. Dopaminergic and adrenergic partial agonist.
Cognition enhancer.

Citations: 2

[View Online »](#)

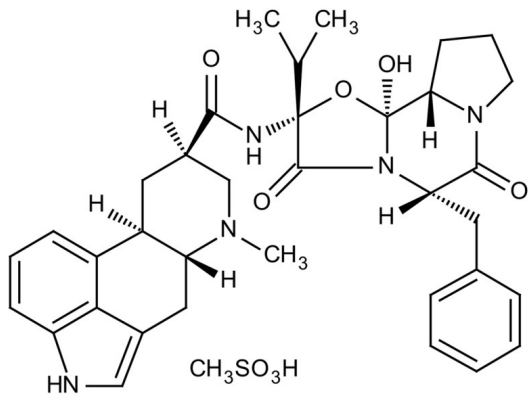
Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| BML-NS108-0200 | 200mg |
|----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 24730-10-7

Couple Target Serotonin receptor

Couple Type Ligand

Formula $C_{35}H_{41}N_5O_5 \cdot CH_3SO_3H$

MW 707.8

Purity $\geq 98\%$

Solubility Soluble in water (10mg/ml), methanol (22mg/ml), or DMSO (25mM).

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)