

# LE300

Dopamine receptor D<sub>1</sub> antagonist

Potent and selective dopamine receptor D<sub>1</sub> antagonist.

Citations: 1

[View Online »](#)

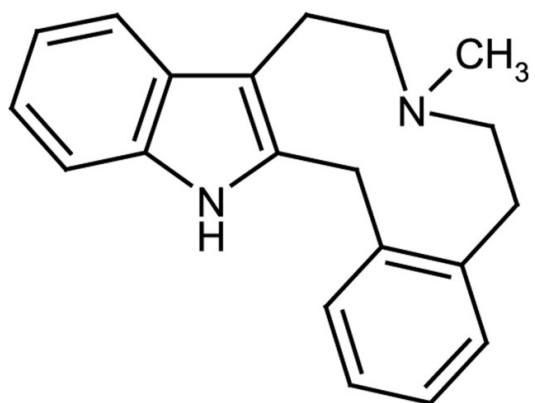
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-550-245-M005 | 5mg |
|------------------|-----|

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                  |
|-------------------|------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Protect from light.                                              |
| Long Term Storage | +4°C                                                             |
| Shipping          | Blue Ice                                                         |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                 |
|------------------|-----------------------------------------------------------------|
| Alternative Name | 7-Methyl-6,7,8,9,14,15-hexahydro-5H-benz[d]indolo[2,3-g]azecine |
| CAS              | 274694-98-3                                                     |
| Couple Target    | Dopamine                                                        |
| Couple Type      | Inhibitor                                                       |
| Formula          | $C_{20}H_{22}N_2$                                               |
| MW               | 290.4                                                           |

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)