

Apomorphine . hydrochloride . hemihydrate

Dopamine agonist

Archetypal dopamine pan-receptor agonist. Displays anti-Parkinsons activity in vivo. Protects against MPTP-induced neurotoxicity in a mouse model. In clinical use for Parkinson's disease.

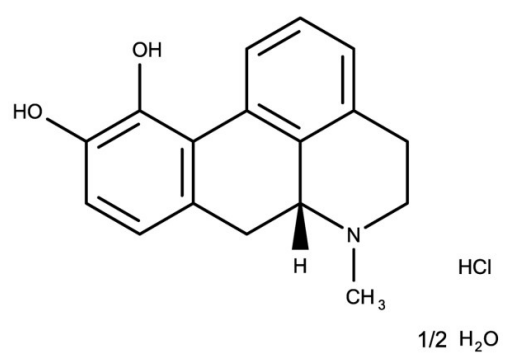
Ordering Information

[Order Online »](#)

ENZ-CHM480-0500	500mg
-----------------	-------

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

Alternative Name (R)-5,6,6a,7-Tetrahydro-6-methyl-4H-dibenzo[de,g]quinolone-10,11-diol hydrochloride

Appearance Grey solid.

CAS 41372-20-7

Couple Target Dopamine receptor

Couple Type Activator

Formula $C_{17}H_{17}NO_2 \cdot HCl \cdot \frac{1}{2} H_2O$

Identity Determined by NMR.

MW 312.78

Purity $\geq 98\%$ (TLC)

Solubility Soluble in water (20 mg/ml).

Last modified: February 27, 2025



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-be@enzolifesciences.com)

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