

8-Hydroxy-DPAT . hydrobromide

5-HT_{1A} agonist

Standard selective 5-HT_{1A} serotonin receptor agonist.

Citations: 3

[View Online »](#)

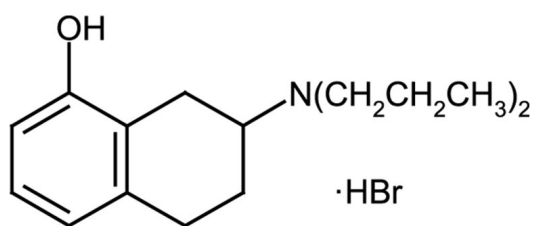
Ordering Information

[Order Online »](#)

ALX-550-195-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	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-8-Hydroxy-2-dipropylaminotetralin . HBr, 8-OH-DPAT . HBr
Appearance	White solid.
CAS	87394-87-4
Couple Target	Serotonin receptor
Couple Type	Inhibitor, Ligand
Formula	C ₁₆ H ₂₅ NO . HBr
MW	328.3
Purity	≥99%
Solubility	Soluble in water or 100% ethanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com