

trans-7-Hydroxy- PIPAT . maleate

Dopamine receptor ligand

A selective, high affinity D₃ dopamine receptor ligand ($K_i=0.99$ nM).

Citations: 1

[View Online »](#)

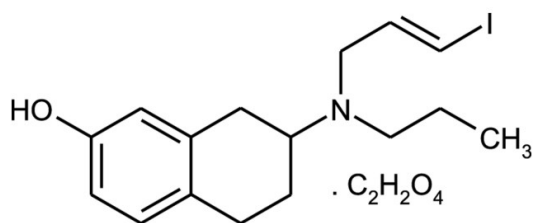
Ordering Information

[Order Online »](#)

BML-D106-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name trans-7-Hydroxy-2-[N-propyl-N(3'-iodo-2'-propenyl)amino]tetralin . maleate

Appearance White solid.

CAS 159559-71-4

Couple Target Dopamine receptor

Couple Type Ligand

MW 487.3

Purity ≥98%

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-fr@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