

Serotonin . hydrochloride

Neurotransmitter

Neurotransmitter. Endogenous 5-HT agonist produced by serotonergic neurons in the CNS and enterochromaffin cells in the gastrointestinal tract.

Citations: 3

[View Online »](#)

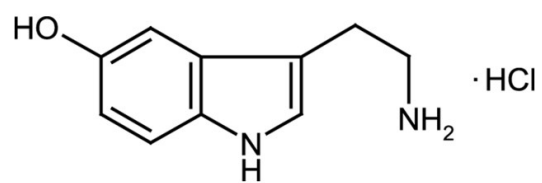
Ordering Information

[Order Online »](#)

ALX-550-328-M050	50mg
ALX-550-328-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Handling Protect from light. Hygroscopic.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-Hydroxytryptamine . HCl

Appearance White to gray crystalline powder.

CAS 153-98-0

Couple Target Serotonin receptor

Couple Type Activator, Ligand

Formula $C_{10}H_{12}N_2O \cdot HCl$

MI 14: 8462

MW 176.2 . 36.5

Purity ≥98% (HPLC)

Solubility Soluble in water, methanol or 100% ethanol.

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-
fr@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)