

Melatonin

Ultra-pure

Endogenous agonist at melatonin MT1 and MT2 receptors. Displays antioxidant and antiapoptotic effects in numerous cell types as well as beneficial effects in various neurodegenerative diseases. It protects neuronal cells from amyloid beta-mediated oxidative damage and increases survival in an Alzheimer's disease transgenic mouse model. Melatonin directly inhibits mitochondrial permeability transition pore ($IC_{50} = 0.8M$) and reduced the loss of neurons in a rat model of transient brain ischemia.

Citations: 21

[View Online »](#)

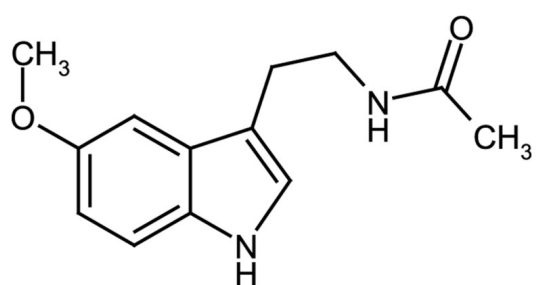
Ordering Information

[Order Online »](#)

BML-NS520-0001	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Acetyl-5-methoxytryptamine
Appearance	White to off-white powder.
CAS	73-31-4
Couple Target	Melatonin receptor, MPTP
Couple Type	Inhibitor, Ligand
Formula	$C_{13}H_{16}N_2O_2$
MI	14: 5816
MW	232.3
Purity	≥98% (HPLC)
RTECS	AC5955000
Solubility	Soluble in dilute aqueous acid, methanol or 100% ethanol.



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)