

# R(-)-Deprenyl . HCl

## Anti-apoptotic and neuroprotective

Deprenyl inhibits apoptosis and displays neuroprotective effects in a variety of *in vitro* and *in vivo* systems. *In vivo* it rescues nigral dopaminergic neurons after systemic MPTP treatment. *In vitro* it rescues PC12 cells from apoptotic cell death induced by trophic withdrawal. The putative target responsible for the anti-apoptotic effects is glyceraldehyde-3-phosphate dehydrogenase. Deprenyl also inhibits monoamine oxidase-B, which is not involved in its anti-apoptotic mechanism. Parkinson's disease therapeutic.

Citations: 4

[View Online »](#)

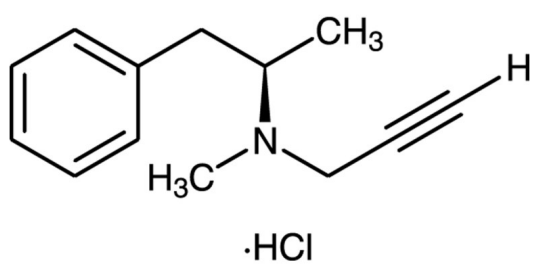
## Ordering Information

[Order Online »](#)

BML-EI240-0100	100mg
----------------	-------

## 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** Selegiline hydrochloride

**Appearance** White solid.

**CAS** 14611-52-0

**Couple Target** GAPDH, Monoamine oxidase

**Couple Type** Inhibitor

**Formula**  $C_{13}H_{17}N \cdot HCl$

**MW** 223.7

**Purity**  $\geq 98\%$  (HPLC)

**Solubility** Soluble in water.

**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)