

β-Synuclein (human), (recombinant)

β-Synuclein co-localizes with α-synuclein in some brain regions but, unlike α-synuclein, it is not found in Lewy bodies of Parkinson's disease. It has been shown to inhibit the self aggregation of α-synuclein.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE258-0500	500μg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Useful for functional studies and drug screening.

Formulation Lyophilized from 10mM Tris-HCl, pH 7.4

MW 14.3 kDa

Source Produced in *E. coli*. full length protein.

UniProt ID Q16143

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-
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)