

[R112A, R113G]Brain derived neurotrophic factor precursor (mouse), (recombinant)

BDNF is a neurotrophic factor expressed by proteolytic cleavage of its precursor, proBDNF. The precursor form was thought to be important for the correct folding, secretion and trafficking of the mature protein.

Trk ligand

Processes such as neuronal process retraction and neuronal apoptosis were thought to be mediated by the binding of BDNF to p75. The finding that proBDNF, and not mature BDNF, is the preferred ligand for p75 has led to a reexamination of the biological roles of the two forms. The amino acid sequence of proBDNF-mut has been changed to prevent the proteolytic cleavage which results in the formation of mature BDNF.

Ordering Information

[Order Online »](#)

BML-SE548-0005	5µg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name pro-BDNF

Application Notes Can be used as growth and differentiation factor for neuronal cultures and for receptor binding studies.

Purity $\geq 99\%$

Source Produced in *E. coli*.

UniProt ID P21237



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)