

# Brain derived neurotrophic factor (human), (recombinant)

Brain derived neurotrophic factor (BDNF) is a member of the NGF family of neurotrophic growth factors and is a prosurvival factor induced by cortical neurons that is necessary for survival of striatal neurons in the brain. It is also required for differentiation of cortical and hippocampal GABAergic interneurons and induces differentiation of dopaminergic neurons and survival of some neuronal cultures. BDNF is expressed as a larger pro form and is cleaved by proteases into its mature form: a 27 kDa dimer formed by two identical subunits.

## Ordering Information

[Order Online »](#)

BML-SE546-0005	5µg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      BDNF

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

Formulation      Lyophilized from a 0.2µm sterile water solution.

Purity      ≥97%

Source      Produced in *E. coli*.

UniProt ID      P23560



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)