

Brain derived neurotrophic factor (human), (recombinant)

Trk ligand
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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com