

# PRODUCT INFORMATION



## BDNF (R129A, R130A mutant; mouse, recombinant)

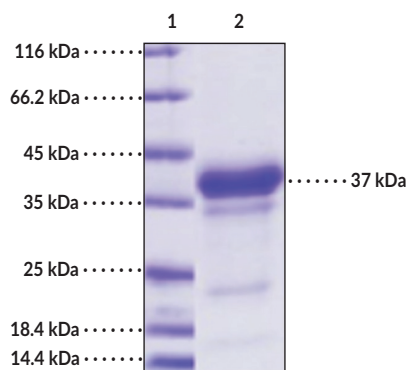
Item No. 42754

### Overview and Properties

**Synonym:** Brain-derived Neurotrophic Factor  
**Source:** Recombinant mouse C-terminal His-tagged BDNF (R129A, R130A mutant) expressed in HEK293 cells  
**Amino Acids:** 19-249  
**Uniprot No.:** P21237  
**Molecular Weight:** 27.4 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥95% estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile PBS, pH 7.4  
**Endotoxin Testing:** <1.0 EU/μg, determined by the LAL endotoxin assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: BDNF

**SDS-PAGE Analysis of BDNF.** This protein has a calculated molecular weight of 27.4 kDa. It has an apparent molecular weight of approximately 37 kDa by SDS-PAGE due to glycosylation.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/12/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

BDNF is a neurotrophin that is involved in synaptic plasticity and the growth, proliferation, and survival of neurons.<sup>1,2</sup> It is synthesized as a precursor protein composed of a signal peptide, a prodomain, and mature BDNF sequence, which is secreted from neurons after cleavage of the proprotein.<sup>1</sup> Incorporation of R129A and R130A mutations into *Bdnf* removes the dibasic amino acid motif that is required for processing of proBdnf to mature Bdnf.<sup>3</sup> *Bdnf* is expressed in several brain regions to varying degrees with the highest levels in the hippocampus and cerebral cortex.<sup>4</sup> BDNF homodimerizes and induces neurotropic signaling through tropomyosin-related kinase B (TrkB) and the p75 neurotrophin receptor (p75<sup>NTR</sup>).<sup>1,5</sup> BDNF increases the proliferation of choriocarcinoma cells *in vitro*.<sup>5</sup> It also induces long-term potentiation (LTP) in rat hippocampus.<sup>6</sup> Knockout of *Bdnf* decreases coordination and balance and sensory neuronal survival in mice.<sup>7</sup> Serum levels of *BDNF* are decreased in patients with depression or schizophrenia and brain expression of *BDNF* mRNA is decreased in patients with Alzheimer's disease.<sup>2</sup> Cayman's BDNF (R129A, R140A mutant; mouse, recombinant) protein consists of 242 amino acids, has a calculated molecular weight of 27.4 kDa, and a predicted N-terminus of Ala19 after signal peptide cleavage. By SDS-PAGE, the apparent molecular mass of the protein is 37 kDa due to glycosylation.

## References

1. Bothwell, M. NGF, BDNF, NT3, and NT4. *Neurotrophic Factors. Handbook of Experimental Pharmacology* Lewin, G. and Carter, B., editors, Springer (2014).
2. Hong, C.-J., Liou, Y.-J., and Tsai, S.-J. Effects of BDNF polymorphisms on brain function and behavior in health and disease. *Brain Res. Bull.* **86(5-6)**, 287-297 (2011).
3. Gaub, P., de Léon, A., Gibon, J., *et al.* HBpF-proBDNF: A new tool for the analysis of pro-brain derived neurotrophic factor receptor signaling and cell biology. *PLoS One* **11(3)**, e0150601
4. Hofer, M., Pagliusi, S.R., Hohn, A., *et al.* Regional distribution of brain-derived neurotrophic factor mRNA in the adult mouse brain. *EMBO J.* **9(8)**, 2459-2464 (1990).
5. Kawamura, N., Kawamura, K., Manabe, M., *et al.* Inhibition of brain-derived neurotrophic factor/tyrosine kinase B signaling suppresses choriocarcinoma cell growth. *Endocrinology* **151(7)**, 3006-3014 (2010).
6. Ying, S.W., Futter, M., Rosenblum, K., *et al.* Brain-derived neurotrophic factor induces long-term potentiation in intact adult hippocampus: Requirement for ERK activation coupled to CREB and upregulation of Arc synthesis. *J. Neurosci.* **22(5)**, 1532-1540 (2002).
7. Ernfors, P., Lee, K.-F., and Jaenisch, R. Mice lacking brain-derived neurotrophic factor develop with sensory deficits. *Nature* **368(6467)**, 147-150 (1994).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM