

GIBCOBRLLIFE  TECHNOLOGIES™

Recombinant Human Brain-Derived Neurotrophic Factor (BDNF)

Cat. No.: 10908-010

Size: 5 µg

Store at -20°C.

Description:

Recombinant Human Brain-Derived Neurotrophic Factor, from *E. coli*, is a 27-kDa-dimer, formed by two identical 119-amino acid subunits, which is held together by strong hydrophobic interactions. It is a potent neurotrophic factor which supports the growth and survivability of nerve and/or glial cells (1-4). For most *in vitro* applications, BDNF exerts its biological activity in the concentration range of 0.1 to 1 ng/ml. Recombinant Human BDNF is provided as a lyophilized protein obtained from a membrane-filtered (0.22 µm) solution of distilled water with no additives.

Directions for Use:

1. Reconstitute in distilled water or aqueous buffers to a concentration of 100 ng/µl. **Note: Do not store in glass;** store in polypropylene vial. Store at -20°C reconstituted for up to six months. Avoid repeated freeze-thaw cycles.
2. The reconstituted solution can be further diluted in distilled water or aqueous buffers to working concentration. **Do not store dilute solutions.**

Note: To avoid adsorptive loss, dilute solutions should be supplemented with Bovine Serum Albumin (BSA) at a concentration of not less than 0.1%.

Quality Control:

- Purity: ≥96% (by SDS-PAGE)
- Endotoxin: ≤ 0.1 ng/µg of BDNF
- Biological Activity: ED₅₀ ≤ 50 ng/ml (determined by dose-dependent induction of choline acetyl transferase activity in primary septal cell cultures)

Doc. Rev.: 031000

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE™ (800) 828-6686.

References:

1. Barde, Y-A, *et al.* (1982) *EMBO J* 1: 549.
2. Martin-Zanca, D. *et al.* (1986) *Nature*. 319: 743.
3. Kalcheim, C. *et al.* (1992) *PNAS* 89: 1661.
4. Snider, W.D. (1994) *Cell* 77: 627.