

PRODUCT DATA SHEET



proBDNF (mouse, recombinant)

CATALOG NO.: SE-547

LOT NO.: Template

DESCRIPTION: proBrain-Derived Neurotrophic Factor (proBDNF) is a 52.4 kDa dimer expressed in *E. coli*.

PURITY: >99% by HPLC.

BIOASSAY: proBDNF was bioassayed for neurotrophic activity in TrkB transfected PC12 cells. Neurite outgrowth was seen from 0.2 nM after 4 days.

SUPPLIED AS: Lyophilized solid. proBDNF was lyophilized from sterile, distilled H₂O.

RECONSTITUTION: proBDNF should be diluted in sterile water or DMEM at a concentration of at least 10 µg/ml.

STORAGE: Lyophilized form: 2-3 weeks at room temperature. 6-9 months at -20° C. After reconstitution: Up to one week at 4° C and 4-6 weeks at -70° C. Avoid repeated freeze/thaw.

REFERENCES:

1. M. Bibel and Y.A. Barde *Genes & Dev.* 2000 **14** 2929
2. S.M. Cahoon-Metzger *et al. Dev. Biol.* 2001 **232** 246
3. C.M. Troy *et al. J. Biol. Chem.* 2002 **277** 34295
4. Z.Y. Chen *et al. J. Neurosci.* 2004 **24** 4401
5. M.F. Egan *et al. Cell* 2003 **112** 257
6. R. Lee *et al. Science* 2001 **294** 1945
7. H.K.Teng *et al. J. Neurosci.* 2005 **25** 5455
8. N.H. Woo *et al. Nat. Neurosci.* 2005 **8** 1069
9. B. Fayard *et al. J. Neurosci.Res.* 2005 **80** 18
10. B. Michalski *et al. Mol. Brain Res.* 2003 **111** 148
11. X.F. Zhou *et al. J. Neurochem.* 2004 **91** 704
12. S.J. Mowla *et al. J. Neurosci.* 1999 **19** 2069
13. S.J. Mowla *et al. J. Biol. Chem.* 2001 **276** 12660
14. T. Matsumoto *et al. Nat. Neurosci.* **2008** **11** 131

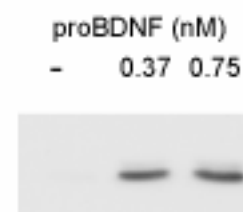


Figure: proBDNF induces ERK1/2 MAPK phosphorylation in neurotrophin receptor tyrosine kinase B (Trk B) transfected HEK-293 cells.

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.