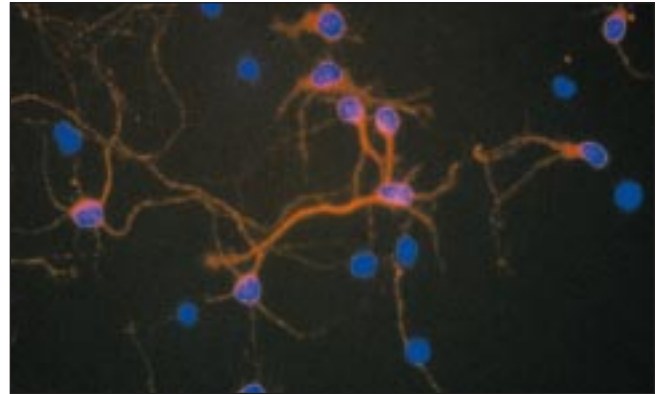


▲ **Anti-NGF mAb, Biotin Conjugate**

Detection of NGF in mouse salivary gland. Details on salivary gland isolation and staining can be found in *Neural Notes* II (3), 11 (1996).

Photograph was kindly provided by Dr. Jean Toma, Montreal Neurological Institute, McGill University, Montreal, Canada.



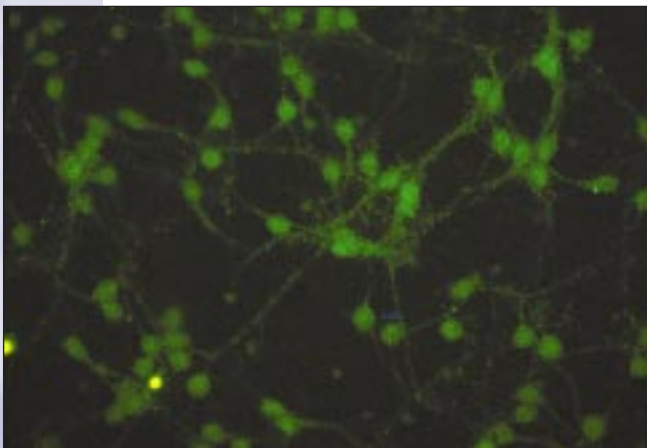
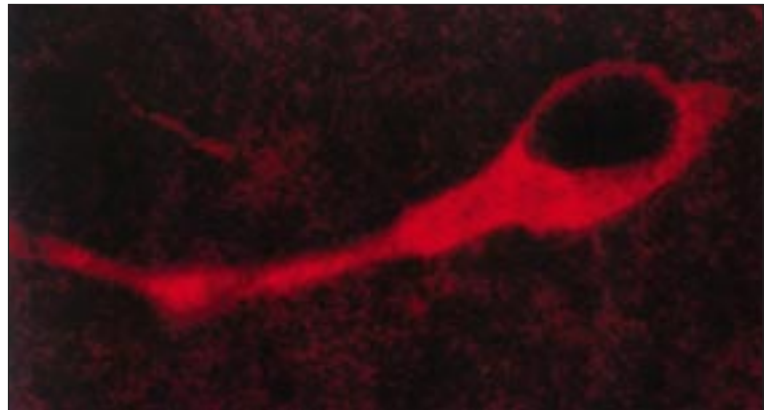
▲ **Anti-Rat CNTF pAb**

Double fluorescence staining of a human fetal CNS cell preparation. Red cells were immunostained with chicken Anti-Rat CNTF pAb and visualized with donkey anti-chicken Cy3 conjugate. Nuclei (blue) are stained with DAPI. Protocol developed and performed at Promega.

Anti-NGF mAb ▶

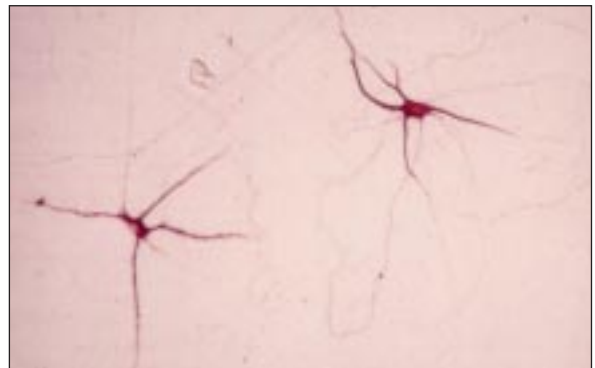
Staining of NGF in a rat basal forebrain cholinergic neuron. Details on neuron isolation and staining can be found on the inside front cover of *Neural Notes* I (4) (1996).

Photomicrograph kindly provided by Dr. Charles Howe, University of California, San Francisco.



▲ **Anti-Human p75 pAb**

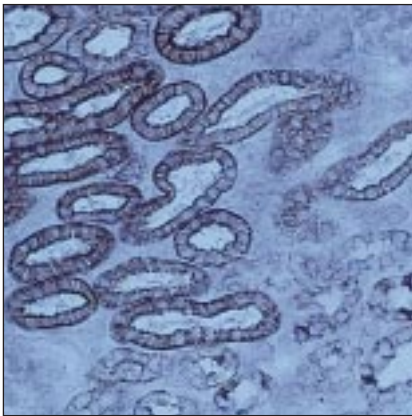
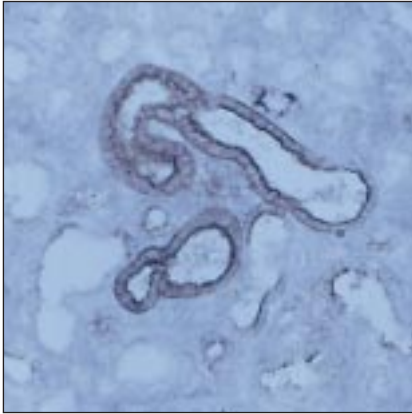
Immunofluorescence staining for p75^{NTR} in a human fetal CNS preparation. A donkey anti-rabbit Cy3 conjugate was used for visualization. Protocol developed and performed at Promega.



▲ **Anti-Human BDNF pAb**

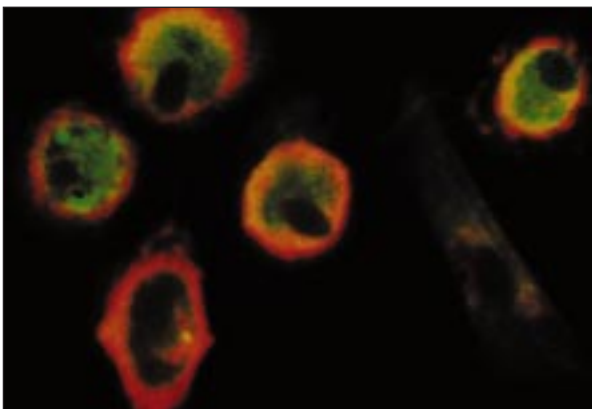
Localization of BDNF in primary cultures of hippocampal neurons. Details on hippocampal neuron isolation, culture and staining can be found in Goodman, L. *et al.* (1996) *Mol. Cell. Neurosci.* 7, 222.

Photomicrograph kindly provided by Dr. Laurie Goodman, Lynx Therapeutics, Hayward, CA. Reprinted by permission of Academic Press.



▲ Anti-TGFβ₁ pAb

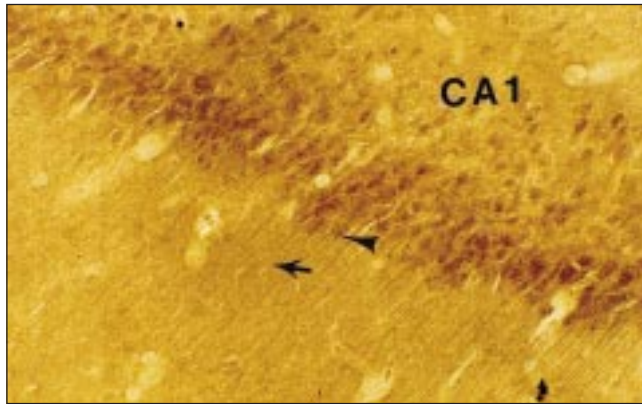
Unfixed adult rat kidney tissue sections stained for TGFβ₁. The image was captured at 200X magnification. Details on tissue isolation and staining can be found in Smith, C. (1995) *Promega Notes* 52, 24.



▲ Anti-Human NT-3 pAb

Immunofluorescent detection of NT-3 in monocyte-derived macrophages and purified human fetal microglia. Microglia cells and macrophages were also reacted against a biotinylated macrophage marker RCA-1 (Ricin Communis Agglutinin 1) that recognizes monocyte derived microglia cells; NT-3 (green) and RCA-1 (red). Details on staining can be found on the inside front cover of *Neural Notes* II (1) (1996).

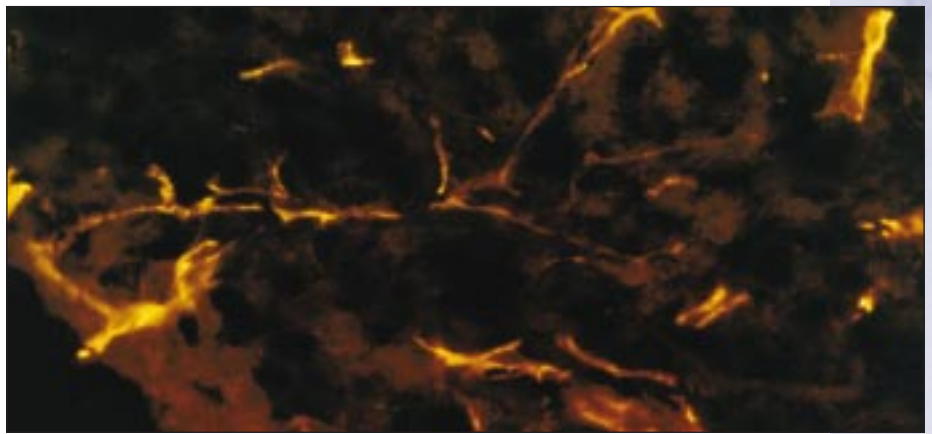
Image kindly provided by Pam Sarnacki, Wanda Wang and Dr. Chris Achim, University of Pittsburgh.



▲ Anti-GBP pAb

Expression of glutamate binding protein in the cell bodies and dendrites of the pyramidal neurons of the hippocampal CA1 region (→ dendrites; ► cell bodies).

Photomicrograph kindly provided by Dr. Eli Michaelis, University of Kansas.



▲ Anti-Pan Trk pAb

Detection of tyrosine kinase (Trk) receptors in rat pineal gland. Details on Anti-Pan Trk pAb can be found in *Neural Notes* II (3), 11 (1996).

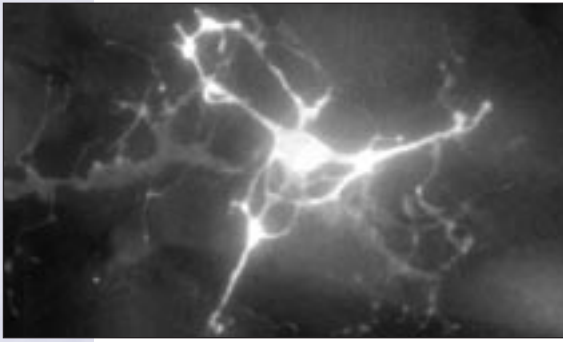
Figure kindly provided by Dr. Jean Toma, Montreal Neurological Institute, McGill University, Montreal, Canada.



▲ Anti-ACTIVE™ MAPK pAb

Detection of the active form of the MAP kinases by immunostaining in PC12 cells. Cells were treated with 200ng/ml of NGF to induce the active form of MAP kinases. Details on Anti-ACTIVE™ MAPK pAb can be found in the following reference *Neural Notes* II (3), 12 (1996).

Photomicrograph kindly provided by Dr. Jean Toma, Montreal Neurological Institute, McGill University, Montreal, Canada.



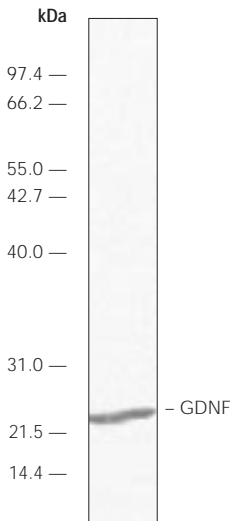
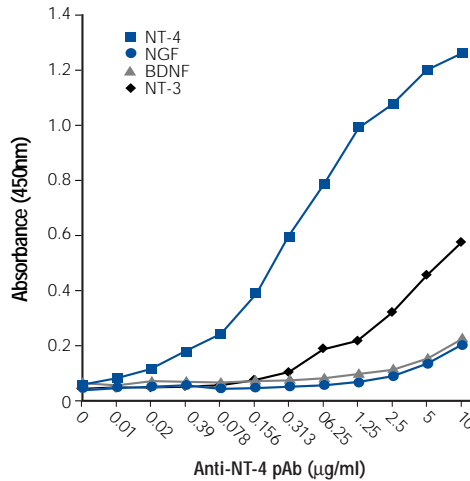
▲ Anti-CNP mAb (Clone 11-5B)

Staining of CNP in rat brain oligodendrocytes. Details on oligodendrocyte isolation, culture conditions and staining can be found in Cammer, W. (1990) *J. Neuroimmunol.* **26**, 173.

Photomicrograph kindly provided by Dr. Wendy Cammer, Albert Einstein College of Medicine, New York.

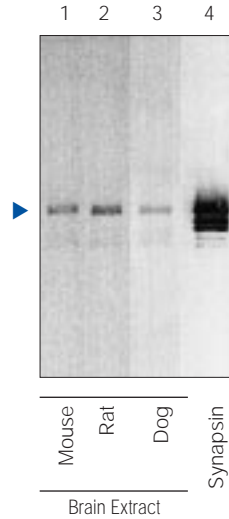
Anti-Human NT-4 pAb

Specificity of Promega's Anti-Human NT-4 pAb demonstrated in an indirect ELISA. Details on the assay procedure can be found in *Neural Notes I* (2), 10 (1995).



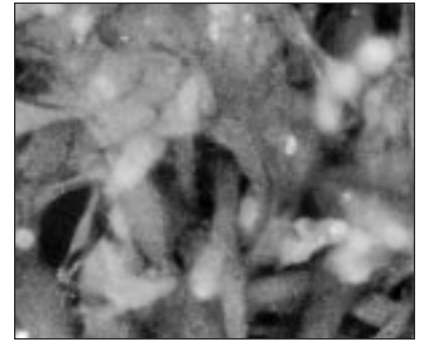
▲ Anti-Human GDNF pAb

Western blot analysis detection of GDNF (100ng). Details on the blotting procedure can be found in *Neural Notes I* (2), 11 (1995).



◀ Anti-Synapsin I N-Terminus mAb

Immunoprecipitation of synapsin from various mammalian brain extracts. Immunoprecipitation also was performed on purified sheep brain synapsin I as shown in lane 4. Details on brain extract preparation and immunoprecipitation can be found in *Neural Notes I* (4), 9 (1996). The arrowhead (▶) indicates the position of synapsin I.



▲ Anti-Human NT-4 pAb

NOBEC-1 cells stained for NT-4. Binding was detected with Anti-Chicken IgY FITC Conjugate (Cat.# G2691). Details on cell culture and staining can be found in *Neural Notes I* (2), 8 (1995).

Antibodies

Product	Size	Cat.#
Anti-Rat CNTF pAb	200µg	G1631
Anti-NGF mAb	100µg	G1131
Anti-NGF mAb, Biotin Conjugate	20µg	G1132
Anti-Human p75 pAb	200µg	G3231
Anti-Human BDNF pAb	200µg	G1641
Anti-TGFβ ₁ pAb	100µg	G1221
Anti-Human NT-3 pAb	200µg	G1651
Anti-Pan Trk pAb	200µg	G1581
Anti-GBP pAb	200µg	G3531
Anti-ACTIVE™ MAPK pAb	15µg	V6671
Anti-CNP mAb (Clone 11-5B)	100µg	G3461
Anti-Synapsin I N-Terminus mAb	1ml	G3351
Anti-Human GDNF pAb	200µg	G2791
Anti-Human NT-4 pAb	200µg	G1621

A complete listing of Promega Antibodies, Neurotrophic Factors and other Growth Factors is available through Promega Online at www.promega.com/neuro/antitab.html