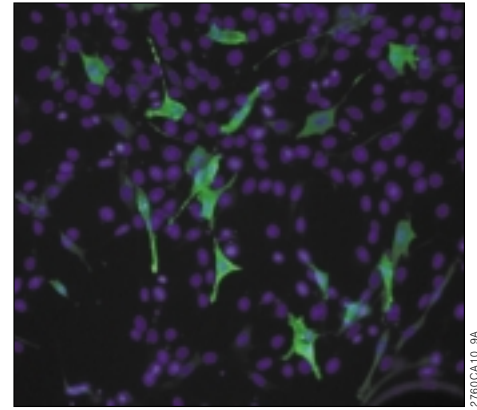


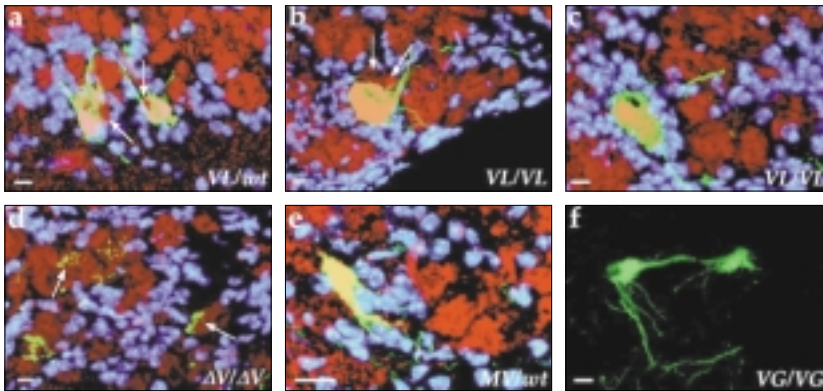
▲ Anti-Cytochrome C mAb

Immunocytochemical labeling for cytochrome C in rat neurosphere (A) and mouse cerebellar granular cell (B) cultures. Cells were stained using Anti-Cytochrome C mAb (clone 6H2.B4, Cat.# G7421). Panel A: E17 rat neurosphere cultures isolated from the subventricular zone striatum and grown in the presence of EGF and serum exhibit typical punctate mitochondrial staining when labeled for cytochrome C (*red*). Dual localization with DAPI (*blue*). Panel B: P8 mouse CGC cultures grown seven days in vitro also show a punctate pattern when immunostained for cytochrome C. Protocols developed and performed at Promega.



▲ Anti-Luciferase pAb

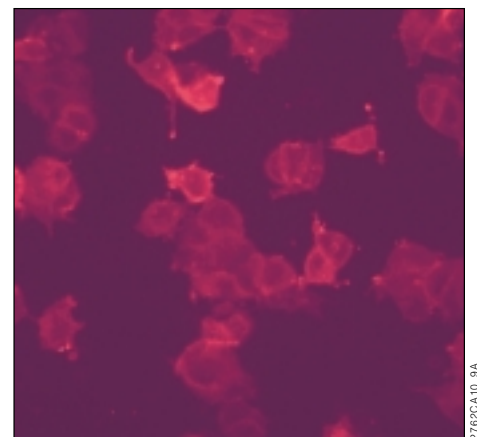
NIH3T3 cells transiently transfected with a luciferase gene. Luciferase-positive cells were detected using Anti-Luciferase pAb (Cat.# G7451). Cells (25,000/well in an ImmunoTech slide chamber) were transiently transfected with a pLuc plasmid using TransFast™ Transfection Reagent^(a) (Cat.# E2431). After 2 days, the cells were fixed using 4% paraformaldehyde, permeabilized with 0.1% Triton® X-100, and blocked with 1% normal donkey serum. Cells were stained with 20µg/ml Anti-Luciferase pAb in PBS for 2 hours followed by 1:200 dilution of donkey anti-goat IgG-FITC (*green*) for 1 hour. Cells were mounted using Vectashield with DAPI (*blue*) and visualized at 200X magnification with a fluorescent microscope and the Spot™ II camera. Protocols developed and performed at Promega.



▲ Anti-β-Galactosidase mAb

Histological analysis of axonal termination in the accessory olfactory bulb (AOB). Sagittal sections through the AOB stained with antibodies against β-galactosidase using Anti-β-Galactosidase mAb (Cat.# Z3781, Z3783; *green*), antibodies against synaptophysin (DAKO, *red*) and DAPI (*blue*). (A) Heterozygous VL mouse; (B, C) Homozygous VL mouse; (D) Homozygous ΔV mouse; (E) Heterozygous MV mouse; (F) Homozygous VG mouse. Details on gene targeting and mutations, as well as culture conditions and immunostaining, may be found in Rodriguez, J., Feinstein, P. and Mombaerts, P. (1999) *Cell* **97**, 199.

Images kindly provided by Dr. Peter Mombaerts, The Rockefeller University, New York. Reprinted by permission of the Cell Press.

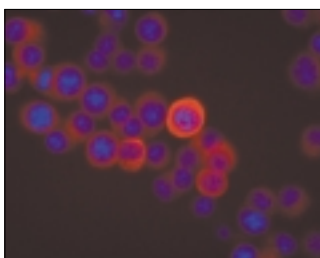


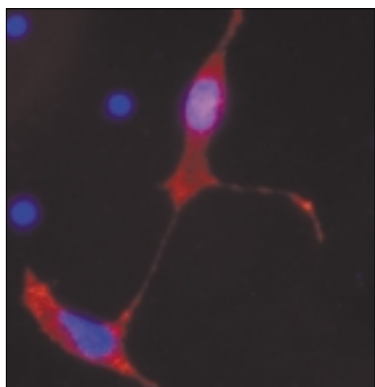
◀ Anti-βIII Tubulin mAb

Undifferentiated PC12 cells immunolabeled for isoform βIII tubulin. βIII tubulin was detected in PC12 cells using 1µg/ml antibody (Cat.# G7121) and visualized with a Cy3™-conjugated secondary antibody (*red*). The DAPI-stained nuclei are visible (*blue*). Oil immersion image, 100X. Protocols developed and performed at Promega.

▲ Anti-Human p75 pAb

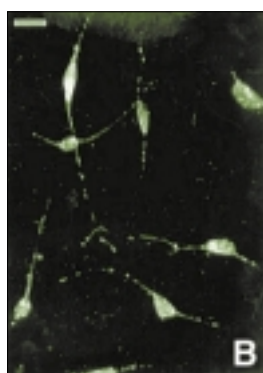
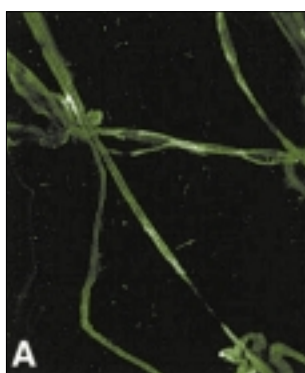
PC12 cells immunolabeled for p75^{NTR}. PC12 cells were incubated with (A) 2µg/ml Anti-Human p75 pAb (Cat.# G3231) or (B) 2µg/ml Anti-Human p75 pAb plus 40µg/ml recombinant GST-p75 and visualized using a donkey anti-rabbit, Cy3™-conjugated secondary antibody (Jackson ImmunoResearch). Protocols developed and performed at Promega.





▲ Immunocytochemical staining of Akt in embryonic rat brain cells
 Embryonic (day 17) rat brain cells were collected and treated with 20ng/ml each of EGF and FGF. Anti-Akt pAb was used at a 1:50 dilution. Positive cells were visualized using a donkey anti-rabbit, Cy3™-conjugated secondary antibody (Jackson ImmunoResearch). Nuclei were stained using DAPI. Protocols developed and performed at Promega.

2766CA10_9A

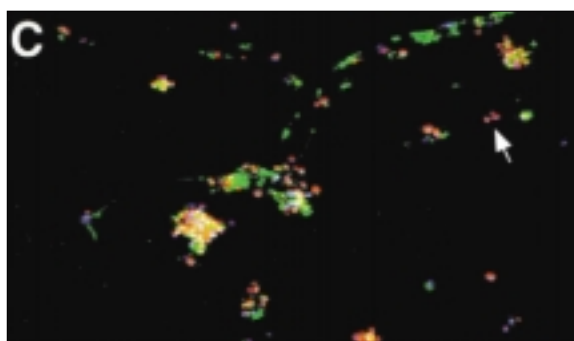
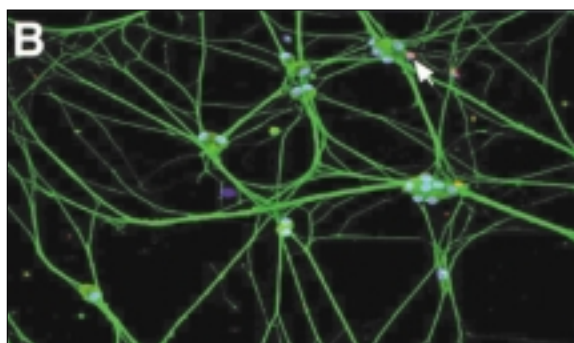
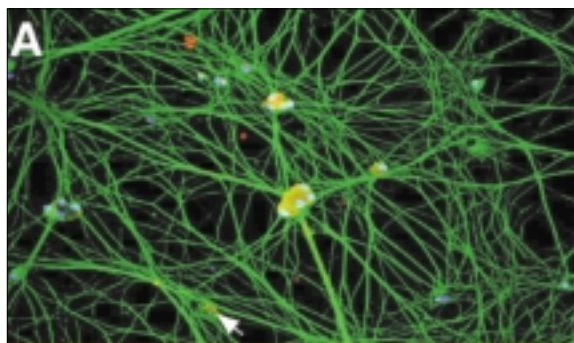


2766CA10_9A

▲ Anti-Human NT-3 pAb

Immunolabeling of NT-3 in teased nerves and cultured Schwann cells. Teased nerves of 4-week-old rats (A) and cultured Schwann cells from 7-day-old (B) rats were immunostained using Anti-Human NT-3 pAb (Cat.# G1651). Details on cell cultures and immunostaining may be found in Meier, C. *et al.* (1999) *J. Neurosci.* **19**, 3847.

Images reprinted by kind permission of Dr. C. Meier and the Society for Neuroscience.



2766CA10_9A

▲ DeadEnd™ Colorimetric Apoptosis Detection System

BDNF decreases neurite outgrowth in cultured neonatal rat sympathetic neurons without decreasing their survival. Digitized micrographs are of postnatal day 1 sympathetic neuron cultures triple-labeled to visualize neurite outgrowth (α-tubulin, green), apoptotic cells (DeadEnd™ System, Cat.# G7130, red/pink) and total number of cells in the culture (Hoechst nuclear stain, blue). Cultures were grown in 50ng/ml NGF for two days and then switched to either (A) 10ng/ml NGF, (B) 10ng/ml NGF plus 100ng/ml BDNF or (C) withdrawn from NGF. Cultures were labeled following an additional two days. Arrows denote nuclei of apoptotic cells. Magnification, 160X. Details on cell cultures and immunostaining may be found in Kohn, J. *et al.* (1999) *J. Neurosci.* **19**, 5393.

Images reprinted by kind permission of Drs. Judi Kohn and Freda Miller, Montréal Neurological Institute, McGill University, Montreal, and the Society for Neuroscience.

Reagents		
Product	Size	Cat. #
Anti-pS473 Akt pAb	40µl	G7441
Anti-Cytochrome C mAb	100µg	G7421
Anti-Luciferase pAb	200µg	G7451
Anti-βIII Tubulin mAb	100µg	G7121
Anti-Human p75 pAb	200µg	G3231
Anti-Human NT-3 pAb	200µg	G1651
Anti-β-Galactosidase mAb	100µg	Z3781
	1mg	Z3783
DeadEnd™ Colorimetric Apoptosis Detection System	40 reactions	G7130
	20 reactions	G7360