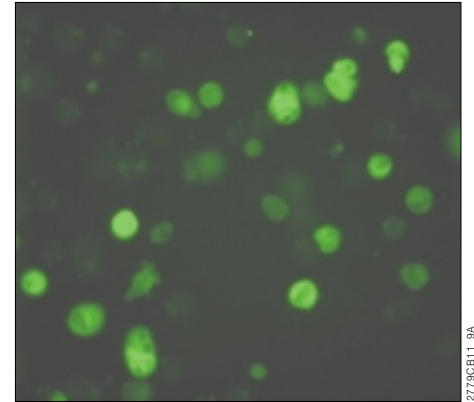


▲ **Anti-Human BDNF pAb**

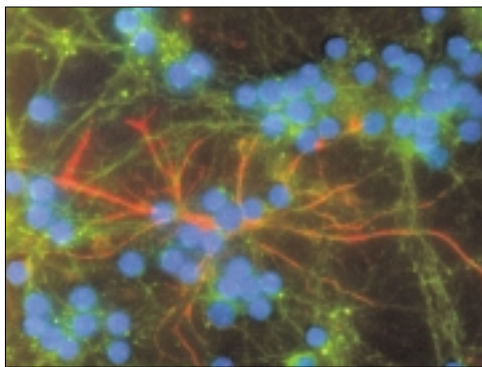
HUVEC cells express factor VIII (von Willebrand factor) and BDNF protein. Low (left) and high (right) magnification of a HUVEC culture stained for BDNF, using Anti-Human BDNF pAb (Cat.# G1641). Additional details on cell cultures and immunostaining may be found in Leventhal, C. *et al.* (1999) *Mol. Cell. Neurosci.* **13**, 450.

Images reprinted by kind permission of Dr. Steven A. Goldman, Cornell University Medical Center, and Academic Press.



▲ **CaspACE™ FITC-VAD-FMK In Situ Marker**

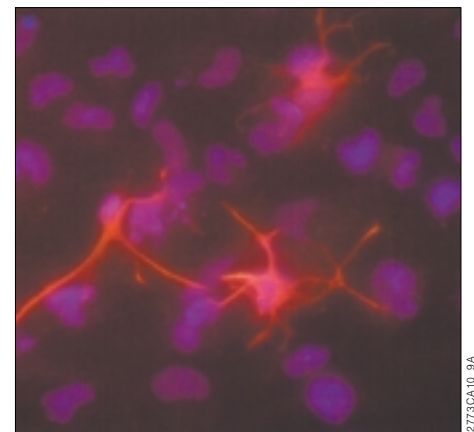
CaspACE™ FITC-VAD-FMK In Situ Marker (Cat.# G7461, G7462) labels anti-Fas mAb-treated Jurkat cells undergoing apoptosis in culture. Protocols developed and performed at Promega.



◀ **Anti-βIII Tubulin mAb**

Immunohistochemical detection of βIII tubulin in mouse cortical granule neurons using Anti-βIII Tubulin mAb (Cat.# G7121). βIII tubulin (green); GFAP (red); DAPI (blue).

Protocols developed and performed at Promega.

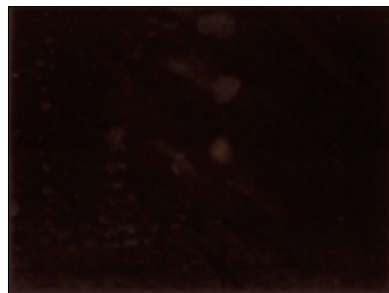
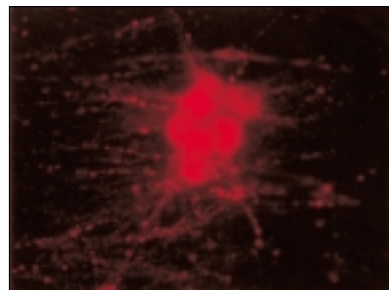
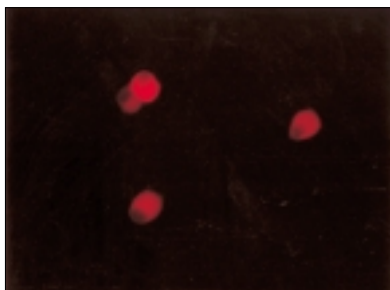


▲ **Anti-GFAP pAb**

Staining of astrocytes by Anti-GFAP pAb. Anti-GFAP pAb (Cat.# G5601) in conjunction with anti-Cy³ conjugate labels astrocytes (red) in mixed population of rat neural progenitor cultures. Nuclei are stained with DAPI (purple). Protocols developed and performed at Promega.

10ng/ml NGF

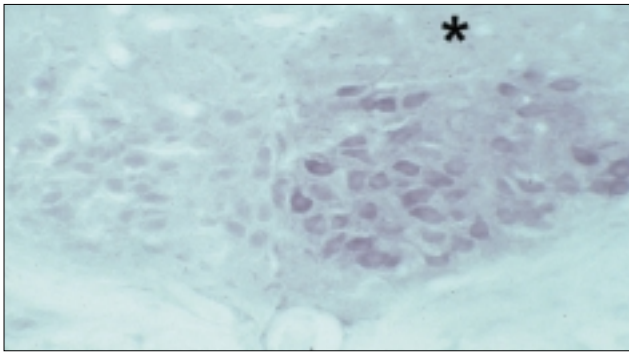
10ng/ml NGF + 50mM KCl



▲ **Anti-ACTIVE® CaM KII pAb**

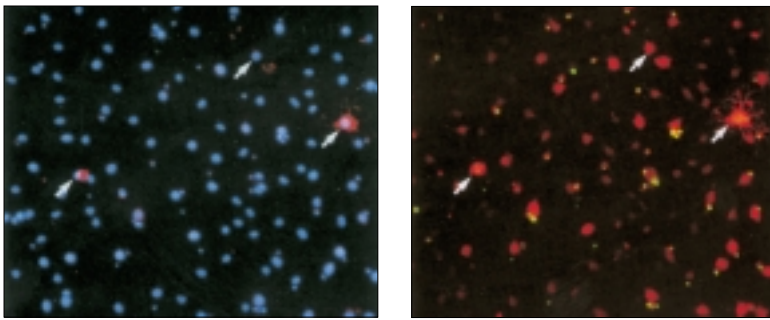
Postmitotic sympathetic neurons were cultured in NGF (10ng/ml; left panels) or NGF plus KCl (50mM; right panels) and treated with KN-62 (10μM; lower panels). KN-62 is a specific pharmacological blocker of CaM KII. Neurons expressing CaM KII were detected using Anti-ACTIVE® CaM KII pAb (Cat.# V1111).

Images kindly provided by Drs. Andrew Vaillant and Freda Miller, Montréal Neurological Institute.



▲ Anti-Human p75 pAb

Immunohistochemical detection of the neurotrophin receptor, p75^{NTR}. Mouse hypoglossal nucleus axotomized on the right side (asterisk) and stained for p75^{NTR}, which is upregulated following axotomy. Anti-Human p75 pAb (Cat.# G3231) was used at a 1:500 dilution. Image kindly provided by Dr. Marie-Pierre Junier, INSERM U421, France.

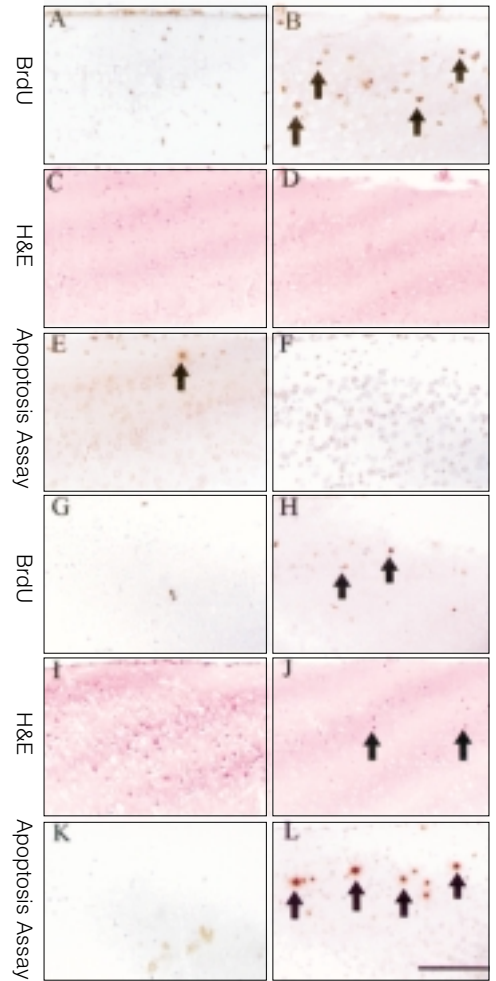


▲ Apoptosis Detection System, Fluorescein

TUNEL analysis of the effect of glial growth factor (GGF) on MBP+ cells. Oligodendrocyte progenitors were differentiated in DM+ culture medium for 5 days and then cultured with or without GGF (200ng/ml) for an additional 18 hours. Immunofluorescence analysis of MBP was combined with TUNEL analysis (Cat.# G3250) of apoptosis cell death. Additional details on cell cultures and immunostaining may be found in Canoll, P.D. *et al.* (1999) *Mol. Cell. Neurosci.* **13**, 79. Images reprinted by kind permission of Dr. James L. Salzer, NYU Medical Center, and Academic Press.

Reagents

Product	Size	Cat.#
CaspACE™ FITC-VAD-FMK In Situ Marker	50µl	G7461
	125µl	G7462
Apoptosis Detection System, Fluorescein	60 reactions	G3250
DeadEnd™ Colorimetric Apoptosis Detection System	40 reactions	G7130
	20 reactions	G7360
Anti-ACTIVE® CaM KII pAb, Rabbit (pT ²⁸⁶)	40µl	V1111
Anti-GFAP pAb	100µg	G5601
Anti-Human p75 pAb	200µg	G3231
Anti-βIII Tubulin mAb	100µg	G7121
Anti-Human BDNF pAb	200µg	G1641



▲ DeadEnd™ Colorimetric Apoptosis Detection System

Cortical sections of wildtype (left) and *Ink4d, Kip1* double-null mice (right) taken at P14 (Panels A–F) and P18 (Panels G–L). At P14, *Ink4d, Kip1* double-null mice show prominent BrdU labeling in layer II of the cortex (Panel B) without apoptosis (Panels D and F), whereas at P18, apoptosis is observed in the cortical layer II (Panels J and L) with less BrdU labeling (Panel H). Bars = 200µm. Details on tissue sectioning procedures, immunohistochemistry and the DeadEnd™ Apoptosis System assay (Cat.# G7130) may be found in Zindy, F. *et al.* (1999) *Proc. Natl. Acad. Sci. USA.* **96**, 13462. Images reprinted by kind permission of Dr. Martine Roussel, St. Jude Children’s Hospital, and the National Academy of Sciences.