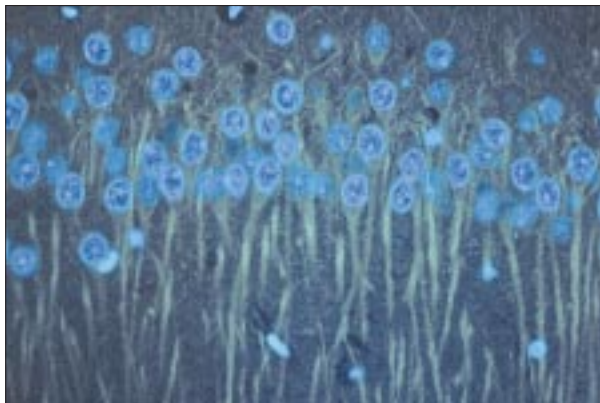
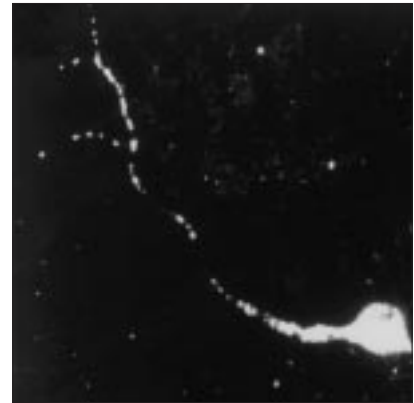


◀ **Anti-βIII Tubulin mAb**

**Immunofluorescence staining of rat cerebellum for βIII tubulin.** **Upper panel:** Frozen rat brain section double-labeled with Anti-βIII Tubulin mAb (anti-mouse Cy3™-conjugated secondary antibody visualized with a Cy3™/Cy5™ filter, red) and DAPI (blue). **Lower panel:** Paraffin-embedded rat brain section; (anti-mouse Cy3™-conjugated secondary antibody visualized with a rhodamine filter, yellow and DAPI, blue). Protocols developed and performed at Promega.

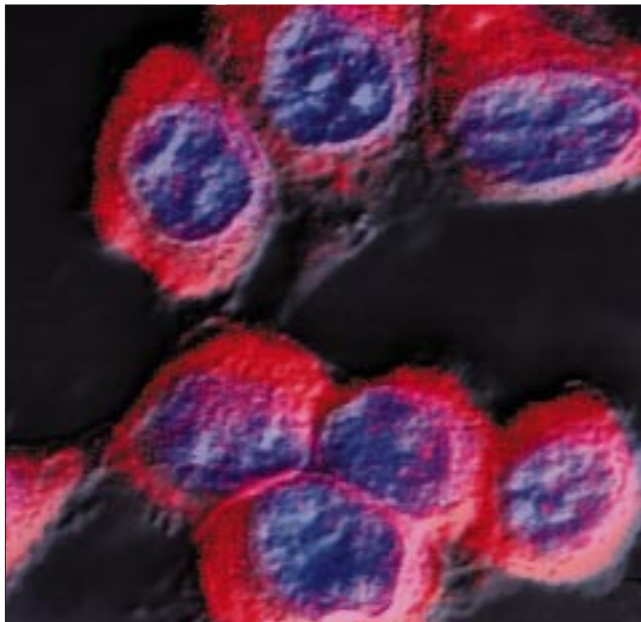
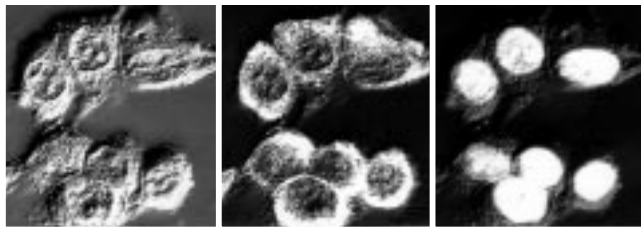
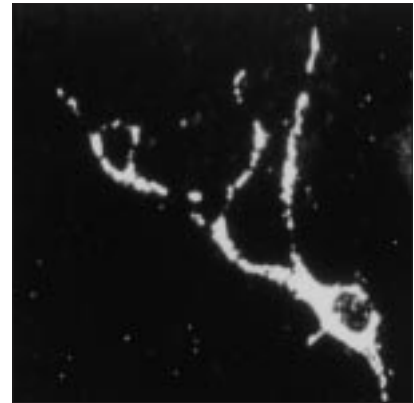


**Anti-Human BDNF pAb** ▶

**Immunocytochemistry of cultured hippocampal neurons incubated in 10mM potassium chloride increases BDNF protein levels in the dendritic compartment.** Anti-BDNF immunostaining of control cultures (**upper panel**); immunostaining following 10mM KCl depolarization (**middle panel**); and

immunostaining after a preincubation with nocodazole for 6 hours followed by KCl treatment and continued incubation in nocodazole (**lower panel**). Details on cell cultures and immunostaining can be found in Tongiorgi, E., Righi, E. and Cattaneo, A. (1997) *J. Neurosci.* **17**, 9492.

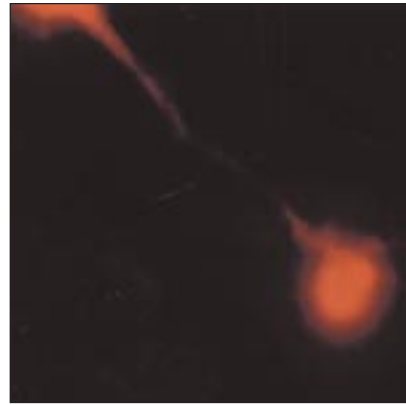
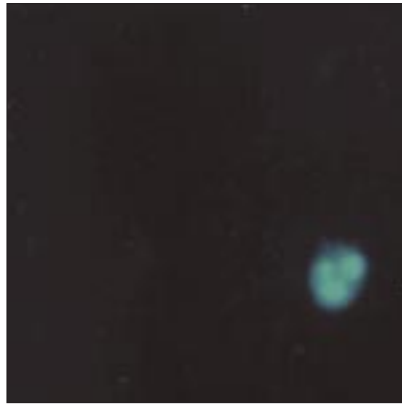
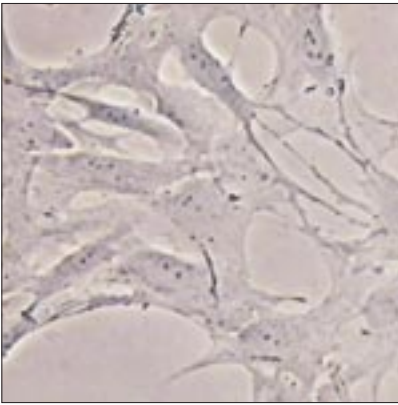
*Images kindly provided by Drs. Enrico Tongiorgi and Antonino Cattaneo, International School for Advanced Studies, Trieste, Italy. Reprinted by permission of The Journal of Neuroscience.*



◀ **Anti-ACTIVE™ MAPK pAb**

**Immunofluorescence analysis of NGF-treated PC12 cells using Anti-ACTIVE™ MAPK pAb.** **Upper panels:** Cells were treated (or not treated, **left panel**) with 50ng/ml NGF for 5 minutes (**middle panel**) or 90 minutes (**right panel**). The upper left panel is a phase contrast image. The images illustrate the translocation of the active form of MAP kinase, for which Anti-ACTIVE™ MAPK pAb is specific, into the nuclei. The **lower image** is a pseudocolor enhancement of the upper middle image; red staining indicates location of active MAP kinase protein.

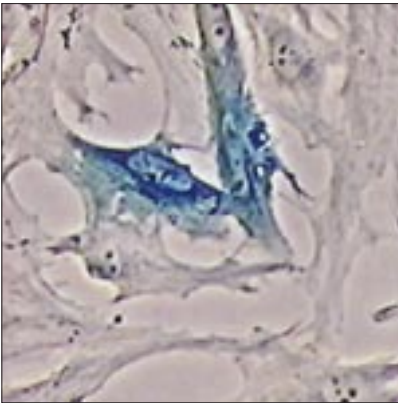
*Images kindly provided by Dr. Philip Lazarovici, Hebrew University, Israel.*



▲ Apoptosis Detection System, Fluorescein

Trophic factor deprivation-induced motor neuron apoptosis. **Left:** Staining of the nucleus of an apoptotic motor neuron cultured for 24 hours without trophic factors. **Right:** Propidium iodide counterstaining of the same motor neuron. Details on the apoptosis assay can be found in Estevez, A.G. *et al.* (1998) *J. Neurosci.* **18**, 923.

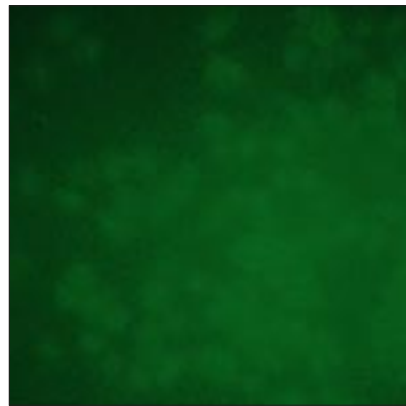
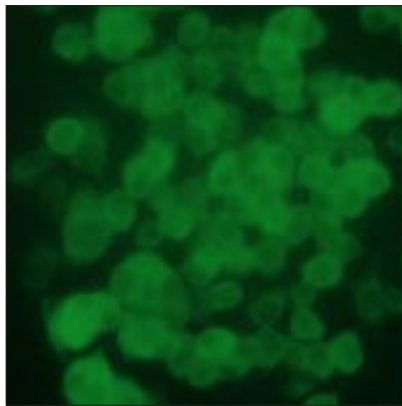
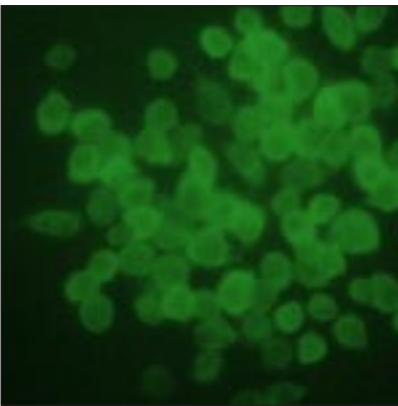
*Images kindly provided by Dr. Alvaro Estevez, University of Alabama at Birmingham. Reprinted by permission of The Journal of Neuroscience.*



◀ TransFast™ Transfection Reagent

Primary culture of mixed glial cells, enriched for astrocytes, from rat cerebellum transfected with TransFast™ Transfection Reagent. A *lacZ* reporter vector was used and positive cells were stained for β-galactosidase expression.

*Photomicrographs kindly provided by Dr. Keith Akama, Northwestern University.*



▲ Anti-ACTIVE™ CaM KII pAb

Immunocytochemical detection of autophosphorylated CaM KII

in PC12 cells. **Left panel:** Cells incubated with antibody only. **Middle panel:** Antibody pre-incubated with a nonphosphorylated CaM KII peptide. **Right Panel:** Antibody pre-incubated with a phosphorylated CaM KII peptide demonstrating specificity for phosphorylated CaM KII. Donkey anti-rabbit, FITC-conjugated secondary antibody (1:500) was used for visualization. Protocols developed and performed at Promega.

Reagents		
Product	Size	Cat.#
Apoptosis Detection System, Fluorescein	60 reactions	G3250
Anti-βIII Tubulin mAb	100µg	G7121
Anti-Human BDNF pAb	200µg	G1641
Anti-ACTIVE™ CaM KII pAb, Rabbit, (pT <sup>286</sup> )	40µg	V1111
Anti-ACTIVE™ MAPK pAb	15µg	V6671
TransFast™ Transfection Reagent*	1.2mg	E2431

\* Patent Pending.

