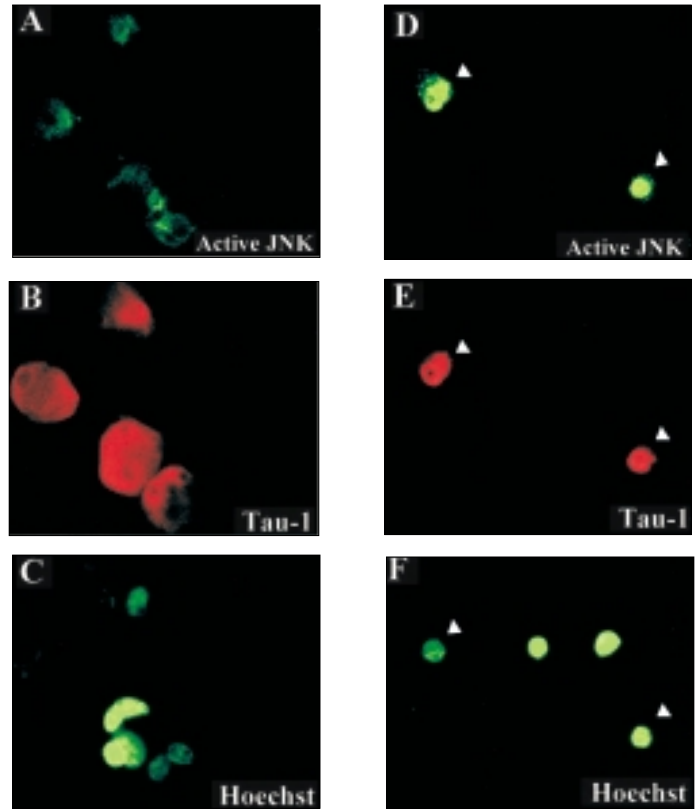


▲ **Anti-Human p75 pAb**

Double-labeled immunofluorescence of neurites extending from an explant plated on Matrigel™. **Panel A:** Immunostaining for β -tubulin III identifies small bundles of neurites extending from the explant. **Panel B:** p75 labeling using Anti-Human p75 pAb (Cat.# G3231) of the same field reveals that olfactory ensheathing cells (OECs) are growing directly below the neurites. Ensheathing cells are spindle-shaped and have aligned in a longitudinal manner to form chains of cells. **Panel C:** Double-label immunofluorescence reveals predominantly yellow-orange olfactory neurites, indicating that they preferentially grow over the surface of OECs rather than the underlying matrix. Details on cell culture and immunostaining may be found in Tisay, K.T. and Key, B. (1999) *J. Neurosci.* **19**, 9890.

Images reprinted by kind permission of Dr. Brian Key, University of Melbourne, and the Society for Neuroscience.

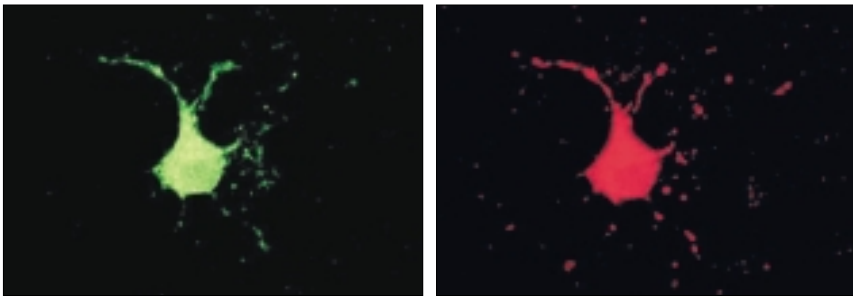


▲ **Anti-ACTIVE® JNK pAb**

Tau-1 staining patterns are altered in apoptotic cells possessing active JNK in the nucleus. Control (+ NGF, Panels A–C) and apoptotic (– NGF, Panels D–F) neuronal PC12 cells deprived of serum for 48 hours were stained with Anti-ACTIVE® JNK pAb (Cat.# V7931, V7932), which recognizes active JNK and was used as a marker of apoptosis, and the tau-1 antibody, which recognizes only an unphosphorylated epitope. Chromatin morphology was assessed using Hoechst stain. Bright nuclear staining of active JNK in cells deprived of serum and NGF indicated apoptosis of neuronal PC12 cells that did not exhibit condensed chromatin (D and F) and that changes in Tau-1 staining patterns were being observed in apoptotic cells (E). Arrowheads point to specific cells (D–F) and are used for orientation purposes between panels. Bar = 5 μ m. Additional details on cultures and immunostaining may be found in Davis, P.K. and Johnson, G.V.W. (1999) *J. Biol. Chem.* **274**, 35686.

Images reprinted by kind permission of Dr. Gail V.W. Johnson, University of Alabama, and The American Society for Biochemistry and Molecular Biology.

2956TA05_0A

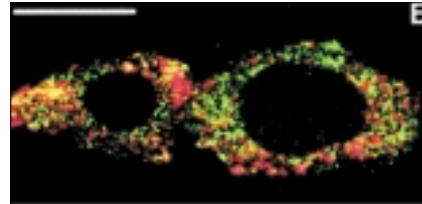
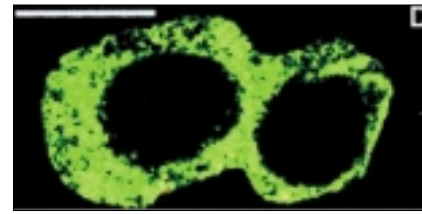


2975CA_06

▲ Anti-CNP mAb

Localization of FasL protein in primary cultures of CNS glial cells by immunohistochemical analysis using polyclonal anti-FasL antibody. Double immunostaining was performed on primary oligodendrocyte cultures using mouse Anti-CNP pAb (Clone 11-5B) (Cat.# G3461; green) and rabbit anti-rat FasL. The mouse antibodies were detected using FITC goat anti-mouse IgG and the rabbit antibody was detected using TRITC goat anti-rabbit IgG. Additional details on cultures and immunostaining may be found in Moalem, G. *et al.* (1999) *FASEB J.* **13**, 1207.

Images reprinted by kind permission of Dr. Michal Schwartz, The Weizmann Institute of Science, and FASEB.



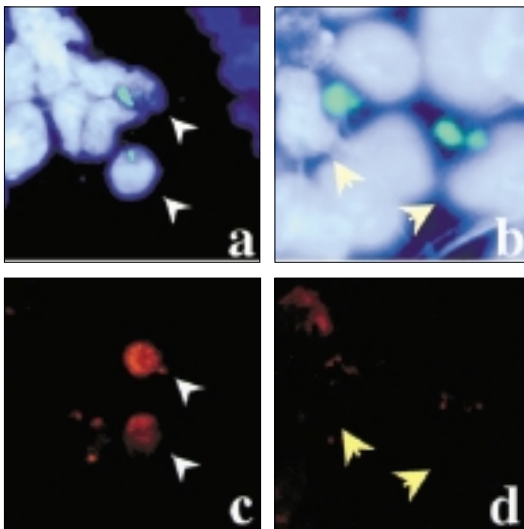
2965TA05_0A

▲ Anti-CNP mAb

Confocal microscopic images of dorsal columns stained to identify oligodendrocytes.

Oligodendrocytes are identified by staining with Anti-CNP pAb (Clone 11-5B) (Cat.# G3461; green). Cytoskeletal damage is demonstrated by marked decrease in spectrin breakdown products (SBP, red) in glutamate-treated (a) versus control (b) slices. Additional details on cultures and immunostaining may be found in Li, S. and Stys, P.K. (2000) *J. Neurosci.* **20**, 1190.

Images reprinted by kind permission of Dr. Peter Stys, University of Ottawa, and the Society for Neuroscience.



2957TA05_0A

▲ DeadEnd™ Colorimetric Apoptosis Detection System

Nuclear aggregate containing cells are preferentially positive in TUNEL assay. TUNEL assay was performed at 48 hours after transfection of tAR97-GFP construct using the DeadEnd™ Colorimetric Apoptosis Detection System (Cat.# G7130, G7360). **Panels A and C:** Arrowheads indicate nuclear aggregate containing cells (Panel A), which were also positive in TUNEL assay (Panel C). **Panels B and D:** Arrows indicate cytoplasmic aggregate containing cells (Panel B), which were negative in TUNEL assay (Panel D) (counterstained with Hoechst 33258). Additional details on cultures and TUNEL assay may be found in Kobayashi, Y. *et al.* (2000) *J. Biol. Chem.* **275**, 8772.

Images reprinted by kind permission of Dr. Gen Sobue, Nagoya University School of Medicine, and the American Society for Biochemistry and Molecular Biology, Inc.

Reagents		
Product	Size	Cat.#
DeadEnd™ Colorimetric Apoptosis Detection System	40 reactions	G7130
	20 reactions	G7360
Anti-ACTIVE® JNK pAb, Rabbit, (pTPpY)	40µl	V7931
	120µl	V7932
Anti-CNP (Clone 11-5B) mAb	100µg	G3461
Anti-Human p75 pAb	200µg	G3231
Anti-βIII Tubulin mAb	100µg	G7121