

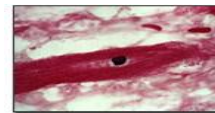
# TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

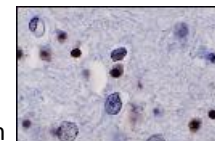
## Specialty TACS® Kits for In Situ Apoptosis Detection

Trevigen's Specialty TACS® Kits were developed to provide researchers with an effective method for identifying apoptotic cells in specific tissue types. These complete kits provide all the reagents required for labeling including permeabilization reagents, labeling and stop buffers, labeling and detection reagents, TACS Nuclease™ reagent, and a protocol detailing incubation times and reagent concentrations.

**CardioTACS™ Kit** – Optimized for identification of apoptotic cells in cardiac samples. The high cellularity of cardiac tissue presents problems in permeabilization, so the kit comes with two permeabilization reagents to provide options. The kit is based on DNA end-labeling using terminal deoxynucleotidyl transferase (TdT) and a modified nucleotide that is subsequently detected using our TACS® Blue Label Detection system.



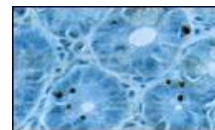
**NeuroTACS™ II Kit** – Optimized to provide rapid and convenient identification of apoptosis in brain tissue or neuronal cells. The kit has been developed to overcome the common difficulties unique to neuronal samples, including the fragile nature of brain tissue sections, high background problems, poor counterstaining with common dyes, and the need to perform dual labeling experiments to detect cell specific antigens in conjunction with apoptotic cells.



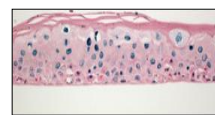
**VasoTACS™ Kit** – Provides an effective method for identifying apoptotic cells in vascular samples, by reducing background and improving labeling in vascular tissue. Similar to the CardioTACS™ Kit, it is based on DNA end-labeling with TdT and a modified nucleotide detected using our TACS® Blue Label Detection System.



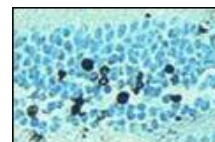
**TumorTACS™ Kit** – Provides rapid and convenient identification of apoptosis in tumors or cancer cells. The kit has a unique TUNEL-based system that preferentially labels the double-stranded DNA found in apoptotic cells. TumorTACS™ contains all the reagents required for permeabilizing, labeling, staining, and counterstaining tumor samples, including TACS-Nuclease™ for preparing positive controls.



**DermaTACS™ Kit** – Optimized for measuring apoptosis in skin samples. The kit is based on DNA end-labeling with TdT and modified nucleotides. Detection of incorporated molecules is achieved using a chromogenic substrate with a horseradish peroxidase detection system. The extracellular scaffolding in skin tissue can make it problematic to permeabilize and high background is a common problem. The DermaTACS™ Kit includes two permeabilization reagents as options for permeabilization along with a detailed protocol with tips for optimal labeling of skin samples.



**TACS.XL® Apoptosis Detection Kit** – For other tissue types than those above, such as the retina tissue shown at right, Trevigen offers the TACS.XL® Kit. The kit's combination of antibody specificity with the signal enhancing properties of biotin: streptavidin result in precise cellular labeling and the highest signal to noise ratio observed in competitive testing.





# TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

## Specialty TACS® Kits for In Situ Apoptosis Detection

### Ordering Information:

Specialty Tissue/Cell Specific TACS® Kits

Description	Size	Catalog #
CardioTACS™ in situ Apoptosis Detection Kit	30 Samples	4827-30-K
NeuroTACS™ II in situ Apoptosis Detection Kit	30 Samples	4823-30-K
VasoTACS™ Kit in situ Apoptosis Detection Kit	30 Samples	4826-30-K
TumorTACS™ Kit in situ Apoptosis Detection Kit	30 Samples	4815-30-K
DermaTACS™ Kit in situ Apoptosis Detection Kit	30 Samples	4829-30-K

### Features and Benefits:

Features	Benefits
Fast assay, streamlined protocol	Requires less than 3 hours to completed
Kits use exclusive TACS® Safe TdT Buffer – sodium cacodylate free	Non-toxic and non-hazardous
Unique buffer system optimized for labeling performance	More consistent labeling
Kits tested on tissue specific samples	Functionally qualified for performance
Only commercially available TUNEL Kits tailored for specific tissue types	Kits include protocol and all reagents necessary for optimal identification of apoptotic cells in specific tissue types

### Related Products:

Description	Size	Catalog #
TACS.XL® Blue Label Kit	30 Samples	4828-30-BK
TACS.XL® DAB Kit	30 Samples	4828-30-DK
TACS.XL® Basic Kit	30 Samples	4828-30-K
TACS.XL® Replenisher Kit	30 Samples	4828-30-R

©2011 Trevigen, TACS and TACS XL are registered trademarks of Trevigen, Inc. CardioTACS, NeuroTACS, VasoTACS, DermaTACS and TACS™ Nuclease are trademarks of Trevigen, Inc.

Trevigen, Inc.  
8405 Helgerman Court  
Gaithersburg, Maryland 20877  
1-800-873-8443