

# TREVIGEN<sup>®</sup> Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

---

## TACS<sup>®</sup> Tissue Control Slides

**Catalog #: 4800-30-40**

**Quantity: 2 slides**

### **Description:**

Trevigen's TACS<sup>®</sup> Tissue Control Slides are ideal for use with any TACS<sup>®</sup> In Situ Apoptosis Detection Kit. The slides provide a sample of immobilized healthy tissue section, allowing users of the TACS<sup>®</sup> In Situ Apoptosis Kits to familiarize themselves with the technique without having to use their own experimental samples. The slides can be used to rule out nonspecific staining and to visualize a small number of apoptotic cells naturally occurring in healthy Thymus. The slides should be treated with TACS-Nuclease<sup>™</sup> to generate the positive control, which will ensure the quality of other components in the TACS<sup>®</sup> Kits.

The paraffin sections (5  $\mu$ m) are from healthy rat thymus tissue, fixed in 3.7% formaldehyde for 24 hours.

**Assay Conditions:** For labeling with TACS<sup>®</sup> In Situ Apoptosis Detection Kits, follow the standard instructions for deparaffinization and rehydration. Permeabilize for 10 minutes at 37°C in a 1 in 50 dilution of proteinase K (cat# 4800-30-01) or in a manner optimized for experimental samples. For a TACS-Nuclease<sup>™</sup> (cat# 4800-30-15) positive control, pretreat the section after permeabilization for 30 minutes at RT with a 1 in 50 dilution of TACS-Nuclease<sup>™</sup> in TACS-Nuclease<sup>™</sup> Buffer (cat# 4800-30-16). TACS<sup>®</sup> Tissue Control Slides can also be used with any of Trevigen's In Situ Apoptosis Detection Kits.

**Storage Conditions:** Store at room temperature with desiccant.

© 2016 Trevigen, Inc. All Rights Reserved. Trevigen, TACS and TACS•XL are registered trademarks and TACS-Nuclease, DermaTACS, NeuroTACS, CardioTACS, TumorTACS, VasoTACS, Apoptosis Grade and Cytonin are trademarks of Trevigen, Inc.

TACS= Trevigen Apoptotic Cell System

E05/09/16v2

---

## TREVIGEN<sup>®</sup>

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

## Related products available from Trevigen:

### Kits

Catalog #	Description	Size
4815-30-K	TumorTACS™ <i>In Situ</i> Apoptosis Detection Kit	30 samples
4823-30-K	NeuroTACS™ <i>In Situ</i> Apoptosis Detection Kit	30 samples
4827-30-K	CardioTACS™ <i>In Situ</i> Apoptosis Detection Kit	30 samples
4829-30-K	DermaTACS™ <i>In Situ</i> Apoptosis Detection Kit	30 samples
4826-30-K	VasoTACS™ <i>In Situ</i> Apoptosis Detection Kit	30 samples
4811-30-K	TACS® 2 TdT Blue Label <i>In Situ</i> Apoptosis Detection Kit	30 samples
4810-30-K	TACS® 2 TdT-DAB <i>In Situ</i> Apoptosis Detection Kit	30 samples
4812-30-K	TACS® 2 TdT Fluorescein Apoptosis Detection Kit	30 samples
4828-30-BK	TACS•XL® Blue Label <i>In Situ</i> Apoptosis Detection Kit	30 samples
4828-30-DK	TACS•XL® DAB <i>In Situ</i> Apoptosis Detection Kit	30 samples

### Antibodies

Catalog #	Description	Size
2291-MC-100	Anti-human-Bcl-2 mAb (clone YTH-8C8)	100 µg
2300-MC-100	Anti-human-Bcl-XL mAb (clone YTH-2H12)	100 µg
2281-MC-100	Anti-human-Bax mAb (clone YTH-6A7)	100 µg
2282-MC-100	Anti-human-Bax mAb (clone YTH-2D2)	100 µg
2305-PC-100	Anti-cleaved human/mouse-Caspase 3 polyclonal	100 µl
4335-MC-100	Anti-PAR polymer mAb (10HA)	100 µl
4336-BPC-100	Anti-PAR polymer polyclonal	100 µl
4338-MC-50	Anti-human/murine-PARP mAb (clone C2-10)	50 µg

### Accessories

Catalog #	Description	Size
4800-30-41	Tissue Control Slides	5 ea.
4864-100	Treated Glass Microscope Slides w/3 sample Hydrophobic Barrier	100 ea.
4867-100	Hydrophobic Coverslips	100 ea.
4862-10	Coverslips 24 x 60 mm, No. 1.5	10 oz.
4865-25	Mounting Medium	25 ml
4800-30-14	Strep-Fluorescein	30 µl
4830-01-3	10X Propidium Iodide	1 ml
4870-500-6	10X PBS	6 x 500 ml
4869-500-6	Apoptosis Grade™ H <sub>2</sub> O	6 x 500 ml
4820-30-13	Blue Counterstain	50 ml

**TACS® Tissue Control Slides (2 ea.)**  
 Catalog #: 4800-30-40  
 Storage: Room Temp.  
**TREVIGEN®**  
 1-800-873-8443