

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## TACS® EpiDerm Control Slides (2 ea.)

**Catalog #:** 4800-30-42

**Contents:** UV irradiated and control EpiDerm sections at 6 µm

**Amount:** 2 slides

**Description:** Trevigen's TACS® EpiDerm Control Slides are ideal for use with the DermaTACS™ In Situ Apoptosis Detection Kit. The sections can be used as a positive control for paraffin embedded tissues and to confirm the quality of other reagents used in combination with TACS® kits. The slides are also useful for the novice who can become familiar with the *in situ* detection of apoptosis without exhausting their own samples. The sample closest to the frosted end of the slide, is untreated EpiDerm and will contain few if any cells with fragmented DNA. The sample farthest from the frosting is EpiDerm at 6 hours following irradiation at 180 mJ/cm<sup>2</sup> and will contain several cells with fragmented DNA. The samples were fixed in 4% formaldehyde for approximately 24 hours and then paraffin embedded. Sections of 6 µm thickness were deposited onto the slides. Typical results obtained using the DermaTACS™ kit are shown below:

### Control EpiDerm

(closest to frosted end)



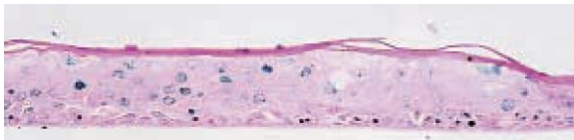
### UV Irradiated EpiDerm

(furthest from frosted end)



### TACS-Nuclease treated

(positive control option)



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

**Assay Conditions:** For labeling with the DermaTACS™ kit, follow the standard instructions for deparaffinization and rehydration. Permeabilize for 30 minutes at 37°C in a 1 in 50 dilution of Proteinase K (cat# 4800-30-01). For a TACS-Nuclease™ (cat# 4800-30-15) positive control, pretreat the section after permeabilization for 10 minutes at 37°C with a 1 in 50 dilution of TACS-Nuclease™ in TACS-Nuclease™ Buffer (cat# 4800-30-16). The EpiDerm Control Slides can also be used with any of Trevigen's In Situ Apoptosis Detection Kits.

© 2010 Trevigen, Inc. All Rights Reserved. Trevigen, TACS and TACS•XL are registered trademarks and TACS-Nuclease, and DermaTACS, NeuroTACS, CardioTACS, TumorTACS, VasoTACS, Apoptosis Grade and Cytonin are trademarks of Trevigen, Inc. EpiDerm is a trademark of MatTek, Inc. TACS= Trevigen Apoptotic Cell System E10/15/10v1

# TREVIGEN®

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 • 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
4411-PC-100	Anti-Phosphorylated Histone-γ-H2AX polyclonal	100 µl
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
2280-MC-100	Anti-mouse-Bax mAb (clone YTH-5B7)	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
6361-PC-100	Anti-human/mouse-PBR 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.
4800-30-21	Cell Culture 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
4825-30-RL	Red Label	30 samples
4878-05-02	Cytonin™ IHC	2 x 5 ml

**TACS® EpiDerm**  
**Control Slides (2 ea.)**  
 Catalog #: 4800-30-42  
 Storage: Room Temp.  
**TREVIGEN®**  
**1-800-873-8443**