

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

TACS® Cell Culture Control Slides (5 ea.)

Catalog #: 4800-30-21

Quantity: 5 slides

Description: The TACS® Cell Culture Control Slides are ideal for use with any of Trevigen's TACS® In Situ Apoptosis Detection Kits. The slides provide a sample of immobilized apoptotic and healthy cells allowing novice users of the TACS® In Situ Apoptosis Kits to familiarize themselves with the technique without having to use their own experimental samples. The slides can be used as a positive control to confirm the quality of other reagents used in combination with the TACS kits.

Area 1 contains apoptotic cells, Area 2 contains untreated cells and Area 3 is empty. Area 1 will show slightly mixed morphology in terms of shape and size and some cell debris may be evident. The counterstaining may be uneven with some cells taking up more counterstain, particularly condensed cells. A majority of cells will be positive for DNA fragmentation indicated by a brown stain (DAB), a blue stain (TACS™ Blue Label) or by bright fluorescence (FITC). Typically, staining is confined to the nucleus although some cells may show spread of positive staining into the cytoplasm caused by small DNA fragments leaking out of the nucleus. Note that apoptotic suspension cells that are stored fixed and dehydrated for long periods of time do not retain good morphology, therefore, interpretation of apoptotic morphology by comparison with these control samples is not recommended. In Area 2 less than 5% cells will be positive for DNA fragmentation. The shape and size of the cells will be similar and should be evenly counterstained.

Sample Preparation: Suspension cells were treated with or without apoptosis inducer, washed with PBS and fixed in 4% formaldehyde. After two washes in PBS, cells were post-fixed in 80% ethanol before being spotted on slides and air dried. After washing in 70% ethanol, the slides were dried at 45°C for 2 hours and stored at 4°C with desiccant.

Assay Conditions: For labeling with TACS® In Situ Apoptosis Detection Kits, the **samples** should be fully rehydrated by passing through a decreasing ethanol series for 5 minutes each (100%, 95%, 70%, 50% ethanol), followed by immersion in PBS for 10 minutes. **For the control slides, permeabilize the sections with Proteinase K for 15 minutes at 37°C. Permeabilized experimental samples and control slide sections should thereafter** be labeled according to the "Instructions For Use" provided with each In Situ Apoptosis Detection kit.

Storage Conditions: Store at 4 °C with desiccant.

© 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
4800-30-42	TACS® EpiDerm Control Slides	2 ea.
4800-30-41	Tissue Control Slides	5 ea.
4800-30-20	Cell Culture Control Slides	2 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 ₂ 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® Cell Culture
Control Slides (5 ea.)**

Catalog #: 4800-30-21

Storage: 4 °C

TREVIGEN®
1-800-873-8443

