

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

TACS[®] Tissue Control Slides (5 ea.)

Catalog #: 4800-30-41

Quantity: 5 slides

Description:

Trevigen's TACS[®] Tissue Control Slides are ideal for use with any TACS[®] In Situ Apoptosis Detection Kit. The slides provide a sample of immobilized healthy tissue section, allowing 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 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[™] to generate the positive control, which will ensure the quality of other components in the TACS[®] Kits.

The paraffin sections (5 μ m) are from healthy rat thymus tissue, fixed in 4% formaldehyde for 24 hours.

Assay Conditions: For labeling with TACS[®] 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[™] (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[™] in TACS-Nuclease[™] Buffer (cat# 4800-30-16). TACS[®] 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.

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
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
4800-30-40	Tissue 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

TACS® Tissue Control Slides (5 ea.)

Catalog #: 4800-30-41

Storage: Room Temp.

TREVIGEN®
1-800-873-8443