

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## TACS•XL® Antibody Module

**Catalog #:** 4828-30-AC

**Size:** 30 samples

**Description:** Apoptosis is characterized by a number of intracellular phenomena such as membrane blebbing, chromatin condensation and nuclear DNA fragmentation. Detection of nuclear DNA fragmentation is a widely accepted method to assay for apoptosis and can be performed *in situ* by incorporating labeled nucleotides onto the free 3' OH ends of DNA fragments using terminal deoxynucleotidyl transferase enzyme (TdT). This type of assay, often referred to as the TUNEL assay, allows monitoring of apoptosis in cell samples or in tissue sections, thus providing histological localization of apoptotic events. TACS•XL® was developed by Trevigen's scientists to improve on background problems that can compromise data interpretation. The TACS•XL® approach utilizes a brominated nucleotide (BrdU) which is more efficiently incorporated by TdT than biotinylated nucleotides at the sites of DNA fragmentation. The incorporated BrdU is detected using a highly specific and sensitive biotinylated anti-BrdU antibody providing superior signal to noise ratios. Our TACS•XL® Antibody Module provides the Anti-BrdU antibody (cat# 4828-30-06) and Strep Diluent (cat# 4828-30-12) for researchers who choose to use their own substrate detection system, or need a reagent for double labeling--using their own antibody(ies) for immunocytochemistry (antigen detection), and our TACS•XL® Antibody for *in situ* detection for apoptosis.

### Antibody solution:

Use 50 µl of diluted antibody per sample. To prepare add:

	2 samples	10 samples	n samples
Anti-BrdU (Cat# 4828-30-06)	2 µl	10 µl	n x 1 µl
Strep-Diluent (Cat# 4828-30-12)	100 µl	500 µl	n x 50 µl

**Assay Conditions:** For labeling instructions for our TACS•XL® In Situ Apoptosis Detection Kits (catalogue numbers: 4828-30-K, 4828-30-BK and 4828-30-DK) please see: <http://www.trevigen.com/apoptosis/tacsXL.php> at [www.trevigen.com](http://www.trevigen.com).

**Storage Conditions:** Store at 4 °C.

## 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](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**TACS™ Cell Culture  
Control Slides (2 ea.)**

**Catalog #: 4828-30-AC**

**Storage: 4 °C**

**TREVIGEN®**

**1-800-873-8443**