

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

TACS•XL™-Replenisher (apoTACS-replenisher)

Catalog #: 4828-30-R

Description: Replenisher reagents for TACS•XL-Basic (Cat# 4828-30-K), TACS•XL-Blue Label (Cat# 4828-30-BK) or TACS•XL-DAB (Cat# 4828-30-DK).

Contents:

Component	Quantity	Storage	Catalog #
Proteinase K Solution	50 µl	-20 °C	4800-30-01
Cytonin	6 ml	4 °C	4876-05-01
B-dNTP Mix	30 µl	-20 °C	4828-30-04
TdT Enzyme	30 µl	-20 °C	4810-30-05
Strep-HRP	30 µl	4 °C	4800-30-06
Anti-BrdU Antibody	30 µl	4 °C	4828-30-06
Strep-Diluent	7.5 ml	4 °C	4828-30-12

Storage: Caution! There are multiple storage temperatures for these products. Please follow appropriate storage temperatures as listed above and on the individual product labels.

Applications: Replenisher reagents for the in situ detection of apoptosis using TACS•XL.

Note: For your convenience, we have included the TACS•XL-DAB as well as the TACS•XL Blue Label complete protocols with your TACS•XL-Replenisher reagents. Please select the appropriate procedure before starting your experiment. If you have any questions, please call our technical service department.

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

**TACS•XL™-Replenisher
(apoTACS-replenisher)**

Catalog #: 4828-30-R

Storage: -20°C

TREVIGEN®
1-800-873-8443