

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

TACS-Nuclease™ and TACS-Nuclease Buffer

Catalog #: 4800-30-N

Contents: TACS-Nuclease **Amount:** 15 µl
 TACS-Nuclease Buffer 1.5 ml

Description: TACS-Nuclease is a highly purified recombinant endonuclease specific for all forms of DNA. TACS-Nuclease has high specific activity, is tolerant to a variety of buffer conditions, and is stable at both 4°C and at -20°C. TACS-Nuclease Buffer has been optimized for use with TACS-Nuclease for the generation of a positive control for DNA fragmentation characteristic of apoptosis. TACS-Nuclease and Buffer are provided as components of Trevigen's TACS® *in situ* apoptosis detection kits:

<u>Kit name</u>	<u>Catalog #</u>
apoTACS™ Blue Label	4828-30-BK
apoTACS DAB	4828-30-DK
CardioTACS™	4827-30-K
NeuroTACS™	4823-30-K
DermaTACS™	4829-30-K
TumorTACS™	4815-30-K
TACS® 2 TdT DAB	4810-30-K
TACS® 2 TdT Blue Label	4811-30-K
TACS® 2 TdT FITC	4812-30-K

Assay Conditions: For use of TACS-Nuclease with any of the TACS *In Situ* Apoptosis Detection Kits listed above, follow the standard instructions provided with the kit for preparation of the samples and for permeabilization. After permeabilization, treat one sample for 10 minutes at 37°C with a 1 in 50 dilution of the TACS-Nuclease prepared in TACS-Nuclease Buffer. Stop the reaction by washing in water or phosphate buffered saline. Note that incubation times up to 60 minutes may be required and empirical testing may be needed for optimal results. After treatment with TACS-Nuclease, proceed with standard labeling, detection and visualization methods as described in the specific kit instructions. Alternatively, include 1 µl of TACS-Nuclease in the standard labeling reaction.

Storage Conditions: May be stored at 4°C or at -20°C. Store at -20°C for maximal stability.

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-Nuclease™ &
TACS-Nuclease™ Buffer**

Cat #: 4800-30-N

Storage: 4 °C or -20 °C
(Manual Defrost Freezer)

TREVIGEN®
1-800-873-8443