

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

TACS® Streptavidin-Horseradish Peroxidase Conjugate (Strep-HRP)

Catalog #: 4800-30-06

Amount: 30 µl

Description: Streptavidin, Mr 60,000, is covalently linked to horseradish peroxidase by a modification of the periodate method with a molar enzyme/protein ratio of 2.5:1. The conjugate binds four molecules of biotin with high affinity ($K_d = 10^{-15} M^{-1}$). The conjugate is provided in liquid form with a proprietary stabilizer added, without preservative.

Source: Streptavidin isolated from *Streptomyces avidinii*

Recommended starting dilutions: TACS® In Situ applications: 1:50 (for DAB based kits) to 1:750 (for TACS Blue Label™ based kits).

Storage Conditions: Store at 4 °C.

Lot Specific Data: see reverse

Related Products available from Trevigen:

Catalog #	Description	Size
4817-60-K	FlowTACS™ Kit	60 samples
4822-96-K	HT TiterTACS™ Assay Kit	96 tests
4830-01-K	TACS® Annexin V FITC Kit	100 samples
4835-01-K	TACS® Annexin V Biotin Kit	100 samples
6300-100-K	DePsipher™ Mitochondrial Potential Assay Kit	100 tests
6305-100-K	MitoShift™ Mitochondrial Potential Assay Kit	100 tests
4815-30-K	TumorTACS™ In Situ Apoptosis Detection Kit	30 samples
4823-30-K	NeuroTACS™ In Situ Apoptosis Detection Kit	30 samples
4829-30-K	DermaTACS™ In Situ Apoptosis Detection Kit	30 samples
4826-30-K	VasoTACS™ In Situ Apoptosis Detection Kit	30 samples
4827-30-K	CardioTACS™ In Situ Apoptosis Detection Kit	30 samples
4828-30-DK	TACS•XL® DAB In Situ Apoptosis Detection Kit	30 samples
4828-30-BK	TACS•XL® Blue Label In Situ Apoptosis Detection Kit	30 samples
4810-30-K	TACS® 2 TdT DAB In Situ Apoptosis Detection Kit	30 samples
4811-30-K	TACS® 2 TdT Blue Label In Situ Apoptosis Detection Kit	30 samples
4812-30-K	TACS® 2 TdT Fluorescein Apoptosis Detection Kit	30 samples
4892-010-K	Cultrex® Calcein-AM Cell Viability Kit	1000 tests
4850-20-ET	TACS® Apoptotic DNA Laddering Kit	20 samples
4684-096-K	HT Colorimetric PARP Apoptosis Assay	96 samples
4685-096-K	HT Chemiluminescent PARP Apoptosis Assay	96 samples

© 2009 Trevigen, Inc. All Rights Reserved. Trevigen, TACS•XL, TACS and Cultrex are registered trademarks, and TACS Blue Label, FlowTACS, TiterTACS, DePsipher, MitoShift, TumorTACS, NeuroTACS, DermaTACS and VasoTACS are trademarks of Trevigen, Inc. TACS = Trevigen Apoptotic Cell System E12/31/09v1

8405 Helgerman Court, Gaithersburg, MD 20877 USA

TREVIGEN®

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

**TACS® Streptavidin-
Horseradish Peroxidase
Conjugate (Strep-HRP)**

Catalog #: 4800-30-06

Storage: 4 °C

TREVIGEN®
1-800-873-8443