



TREVIGEN®

DNA Laddering Kits for Apoptosis Research

The TACS Apoptotic DNA Laddering Kits are used to detect and estimate the level of internucleosomal DNA fragmentation that occurs during apoptosis. Kit selection is dependent upon the degree of apoptosis, number of cells and staining/ detection preference. The evidence of DNA laddering supports other experimental data derived from morphological identification methods. Each kit contains all reagents necessary to isolate, label and detect DNA.

Ordering Information

Product Description	Size	Catalog #
Ethidium Bromide DNA Laddering Kit	20 Samples	4850-20-ET
Isotopic DNA Laddering Kit	20 Samples	4850-20-K
Chemiluminescent DNA Laddering Kit	20 Samples	4855-20-K
Colorimetric DNA Laddering Kit	20 Samples	4857-20-K

Kit Selection Criteria

Product Description	Sensitivity	% Apoptosis	# Cells	Laboratory Requirements
Ethidium Bromide DNA Laddering Kit	Low	20%	1 x 10 ⁶	<ul style="list-style-type: none"> • Photography equipment • EtBr
Isotopic DNA Laddering Kit	Best	1%	1 x 10 ⁵	<ul style="list-style-type: none"> • Southern blotting • Radioisotopes • Darkroom
Chemiluminescent DNA Laddering Kit	Better	2%	1 x 10 ⁵	<ul style="list-style-type: none"> • Southern blotting • Darkroom
Colorimetric DNA Laddering Kit	Good	10%	1 x 10 ⁵	<ul style="list-style-type: none"> • Southern blotting

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS



TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

DNA Laddering Kits for Apoptosis Research

Related Products

TACS•XL® embodies a new approach for the in situ detection of apoptosis. The TACS•XL kit is based on incorporation of bromodeoxyuridine (BrdU) at the 3' OH ends of the DNA fragments that are formed during apoptosis. The incorporation of BrdUTP by TdT is more efficient than either biotinylated or digoxigenin labeled nucleotides used in other TUNEL-based assays. The detection system utilizes a biotin conjugated anti-BrdU antibody and streptavidin-horseradish peroxidase. The combination of antibody specificity with the signal enhancing properties of biotin:streptavidin results in precise cellular labeling and the highest signal to noise ratio observed in competitive testing.

Product Description	Size	Catalog #
TACS.XL® Blue Label Kit	30 Samples	4828-30-BK
TACS.XL® DAB Kit	30 Samples	4828-30-DK
TACS.XL® Basic Kit	30 Samples	4828-30-K
TACS.XL® Replenisher Kit	30 Samples	4828-30-R

The TACS™ 2 TdT Kits utilize Trevigen's unique cation optimization system to enhance labeling within particular tissues. The TACS™ 2 TdT Kits employ a proprietary labeling buffer that contains no toxic components (e.g. sodium cacodylate). A highly purified form of the TdT enzyme is included in the kits for the enzymatic incorporation of biotinylated nucleotides. Biotin labeling is achieved using streptavidin-horseradish peroxidase, and colorimetric substrates diaminobenzidine (DAB) or TACS™ Blue Label™. For fluorescent detection, a fluorescein conjugate of streptavidin is used and visualized by epifluorescence microscopy. Additional fluorophores are now available separately from Trevigen.

Product Description	Size	Catalog #
TACS 2 TdT-Core Kit	30 Samples	4810-30-CK
TACS 2 TdT-Fluorescein Kit	30 Samples	4812-30-K
TACS 2 TdT-Blue Label Kit	30 Samples	4811-30-K
TACS 2 TdT-DAB Kit	30 Samples	4810-30-K

TACS.XL is a registered trademark of Trevigen, Inc.

Copyright 2010