

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

T4 Endonuclease V FLARE[™] Module

Catalog #: 4055-100-FM

Contents:

<u>Component</u>	<u>Catalog #</u>	<u>Amount</u>	<u>Storage</u>
T4-Endo V	4055-100-01	100,000 units	4 °C
25X FLARE [™] Buffer 1	3950-040-01	40 ml	Room temp.
REC [™] Dilution Buffer	3950-010-03	10 ml	-20 °C
100X BSA Additive	3950-100-04	100 µl	-20 °C

Description: Reagents for analysis of DNA damage in single cells using the Trevigen CometAssay[®], FLARE[™] Slides, and T4 endonuclease V.

Assay & Analysis: Refer to the T4 Endonuclease V FLARE[™] Assay Kit's (cat#: 4055-100-FK) Instructions for Use for further details.

Note: CometSlides and FLARE Slides are interchangeable in the FLARE Assay. The difference is in the sample number; the CometSlide[™] holds 2 samples whereas the FLARE Slide holds 3.

Lot Specific Data: Contact Trevigen for lot specific activity of T4 endonuclease V.

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

Related Products:

FLARE™ Assay Kits:

Catalog #	Description	Damage Recognized	Size
4040-100-FK	Fpg Kit	8-oxoguanine, DNA containing	75 samples
4040-100-FM		formamidopyrimidine moieties	100 samples
4055-100-FK	T4 Endo V Kit	Cis-syn isomers of cyclo-	75 samples
		butane pyrimidine dimers	100 samples
4065-100-FK	cv-PDG Kit	Cis-syn and trans-syn isomers of	75 samples
4065-100-FM		cyclobutane pyrimidine dimers	100 samples
4130-100-FK	hOGG1 Kit	8-oxoguanine, DNA containing	75 samples
4130-100-FM		formamidopyrimidine moieties	100 samples
4100-100-FK	UVDE Kit	Cyclobutane pyrimidine	75 samples
4100-100-FM		dimers, (6-4) photoproducts	100 samples
4045-010-FK		Thymine Glycol, 5,6-dihydro-	75 samples
4045-010-FM	III Kit	thymine, urea, 5-hydroxy-6-	100 samples
		hydrothymine, 5,6-dihydro- uracil, alloxan, 5-hydroxy-6- hydrouracil, uracil glycol, 5- hydroxy-5-methylhydantoin, 5-hydroxycytosine, 5-hydroxy uracil, methyltaronyl urea, thymine ring saturated or fragmentation product	

CometAssay® Kits:

Catalog #	Description	Size
4250-050-ESK	CometAssay® Starter Kit	each
4250-050-ES	CometAssay® Electrophoresis System	1 system
4250-050-K	CometAssay®	50 samples
4252-040-K	CometAssay® Higher Throughput Kit	40 samples
4254-200-K	CometAssay® Silver Staining Kit	200 samples
4253-096-K	CometAssay® Kit 96 Wells	96 samples

**T4-Endo V FLARE™
Module**

Catalog #: 4055-100-FM
Storage: 4 °C, -20 °C &
Room Temp.

TREVIGEN®
1-800-873-8443