

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

FLARE[™] Slides

Catalog #: 3950-075-02

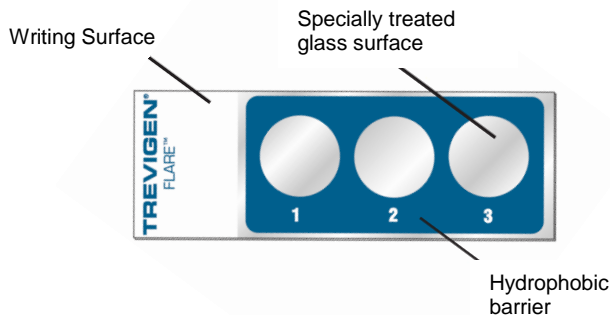
Size: 25 slides

Description: Trevigen's FLARE[™] Slide is a component of the FLARE[™] Kits used for the rapid analysis of DNA adducts revealed by using the comet assay in conjunction with DNA repair enzymes. The comet assay or single cell gel electrophoresis assay permits the quantification of DNA damage in individual cells based on the migration of DNA structures in an electric field. As a result of the migration, comets are formed and the amount of DNA in the comet tail as well as the distance of migration, can be a measure of damage. Trevigen's FLARE[™] Slide provides a convenient, simple and reliable sample preparation, processing, analysis and storage format for the comet and FLARE assay.

Physical State: Sample-ready pretreated glass slides.

Storage: Store at room temperature with desiccant.

Applications: For use with Trevigen's CometAssay[®] and FLARE[™] assay reagents and protocols.



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 Available From Trevigen

FLARE™ Assay Kits:

Catalog #	Description	Damage Recognized	Size
4040-100-FK	Fpg Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4040-100-FM			100 samples
4055-100-FK	T4-PDG Kit	Cis-syn isomers of cyclobutane pyrimidine dimers	75 samples
4055-100-FM			100 samples
4065-100-FK	cv-PDG Kit	Cis-syn and trans-syn isomers of cyclobutane pyrimidine dimers	75 samples
4065-100-FM			100 samples
4100-100-FK	UVDE Kit	Cyclobutane pyrimidine dimers, (6-4) photoproducts	75 samples
4100-100-FM			100 samples
4045-01K-FK	Endonuclease III Kit	Thymine Glycol, 5,6-dihydrothymine, urea, 5-hydroxy-6-hydrothymine, 5,6-dihydro-uracil, alloxan, 5-hydroxy-6-hydro-uracil, uracil glycol, 5-hydroxy-5-methylhy-dantoin, 5-hydroxycytosine,5-hydroxy-uracil, methyl-taronylurea, thymine ring saturated or fragmentation product	75 samples
4045-01K-FM			100 samples
4130-100-FK	hOGG1 Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4130-100-FM			100 samples

CometAssay® Kits:

Catalog #	Description	Size
4250-050-ESK	CometAssay® Starter Kit	each
4250-040-K	CometAssay®	40 samples

Accessories:

Catalog #	Description	Size
4250-050-03	CometSlide™ (2 well)	25 slides
4250-200-03	CometSlide™ (2 well)	100 slides
4252-200-01	CometAssay® HT Slide (20 well)	10 slides
4253-960-03	96 Well CometSlide™	10 slides
4256-010-CC	CometAssay® Control Cells	10 assays
4250-050-ES	CometAssay® Electrophoresis System	each

FLARE™ Slides
 Catalog #: 3950-075-02
 Storage: Room
 Temperature

TREVIGEN®
 1-800-873-8443