

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

CometSlide™

Catalog #: 4250-200-03

Size: 100 slides

Description: Trevigen's CometSlide™ is a component of the CometAssay® Kit (cat# 4250-050-K) used for the rapid analysis of DNA fragmentation associated with DNA damage. 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. Evaluation of the DNA "comet" tail shape and migration pattern allows for assessment of DNA damage, expressed as tail length, percent DNA in the tail and tail moment. Trevigen's CometSlide™ provides a convenient, simple and reliable sample preparation, processing, analysis and storage format for the comet 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.

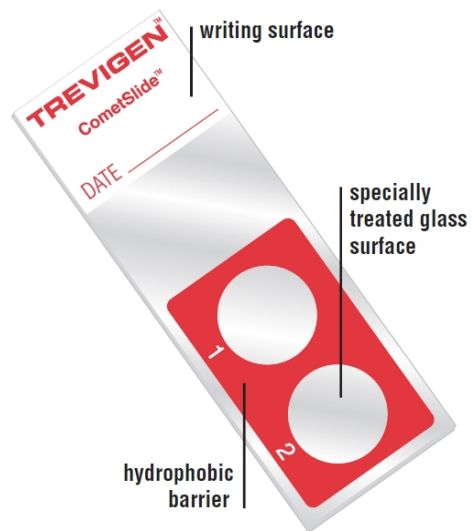


Figure 1. CometSlide™ features

CometAssay® Kits:

Catalog #	Description	Size
4250-050-ESK	CometAssay® Starter Kit	each
4250-050-ES	CometAssay® ES	each
4250-050-K	CometAssay®	50 samples
4251-050-K	CometAssay® Silver Kit	50 samples
4254-200-K	CometAssay® Silver Staining Kit	200 samples
4252-040-K	CometAssay® Higher Throughput Kit	40 samples
4253-096-K	CometAssay® Kit 96 Wells	96 samples

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
4130-100-FK	hOGG1 Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4130-100-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-hydroxy-mine, 5,6-dihydro-uracil, alloxan, 5-hydroxy-6-hydroxy-uracil, uracil glycol, 5-hydroxy-5-methylthio-dantoin, 5-hydroxycytosine, 5-hydroxy-uracil, methyl-taronylurea, thymine ring saturated or fragmentation product	75 samples
4045-01K-FM			100 samples

Accessories:

Catalog #	Description	Size
4256-010-CC	CometAssay® Control Cells (alkaline assay)	1 set
4257-010-NC	Neutral CometAssay® Control Cells	1 set
4252-200-01	CometSlide™ HT (20 well)	10 slides
4253-960-03	96 Well CometSlide™	1 slide
3950-300-02	FLARE™ Slides (3 well)	100 slides

©2011 Trevigen, Inc. All rights reserved. Trevigen and CometAssay are registered trademarks and CometSlide and FLARE are trademarks of Trevigen, Inc. E2/9/11v1

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@travigen.com • www.trevigen.com

CometSlide™
100 Slides

Catalog #: 4250-200-03
 Store at Room Temperature
TREVIGEN®
 1-800-873-8443