

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

CometSlide™ Rack System

Catalog #: 4252-040-02

Description: The CometSlide™ Rack System was designed to accommodate Trevigen's 2 well, 3 well and 20 well CometSlide™ for use in the CometAssay® and FLARE™ assays.

Contents: Each CometSlide™ Rack System is composed of a durable plastic rack that holds 20 slides, and a translucent plastic reservoir with lid.

Instructions: For CometSlide™ (1" x 3"), add 200 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

For the CometSlide™ HT (2" x 3"), add 350 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

Product Care: A mild laboratory detergent may be used, if desired, followed by thorough rinsing with deionized water. The system may be sprayed down with 70% Ethanol for decontamination. Rinse thoroughly with deionized water before each use. **Do not Autoclave! Exposure to ultraviolet light for extended periods of time may compromise integrity of product. Not compatible with acetone or xylenes.**

Storage: Store at Room Temperature.

Related Products Available from Trevigen:

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-tartronylurea, 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
4250-050-03	CometSlide™ (2 well)	25 slides
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

©2010 Trevigen, Inc. All rights reserved. Trevigen and CometAssay are registered trademarks and CometSlide and FLARE are trademarks of Trevigen, Inc. E10/15/10v1

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™ Rack System
Catalog #: 4252-040-02
Store at Room Temperature
TREVIGEN®
1-800-873-8443