

# TREVIGEN<sup>®</sup> Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

---

## CometSlide<sup>™</sup> Rack System

**Catalog #:** 4252-040-02

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

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

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

For the CometSlide<sup>™</sup> 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.

©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<sup>®</sup>

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](mailto:info@travigen.com) • [www.trevigen.com](http://www.trevigen.com)

**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-hydrothy-mine, 5,6-dihydro-uracil, alloxan, 5-hydroxy-6-hydrouracil, uracil glycol, 5-hydroxy-5-methylhy-dantoin, 5-hydroxycytosine,5-hydroxy-uracil, methyl-tartonylurea, 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 |

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