

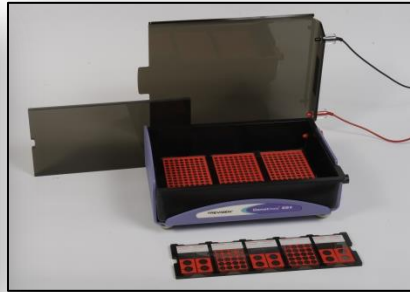
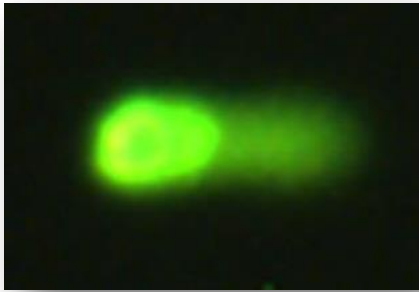


Trevigen®

Better Products. Better Results.

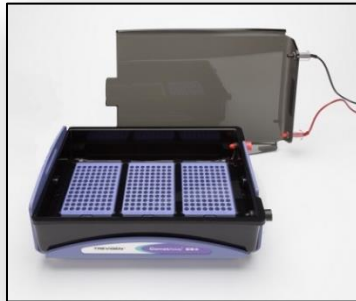
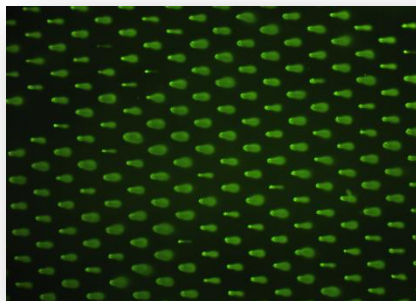
CometAssay® System

Single cell gel electrophoresis or Comet Assay is a technique for quantitating DNA damage and repair from in vivo and in vitro samples of eukaryotic cells. It is the only technique that directly measures DNA damage in individual cells and as a result has gained importance in the fields of genetic toxicology and human biomonitoring. Trevigen offers the complete, optimized system: slides, reagent kits, control cells, electrophoresis unit and analysis software.



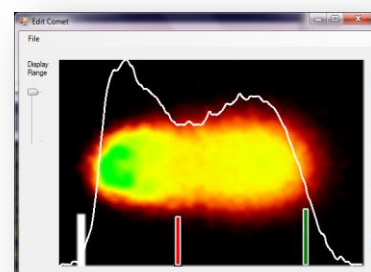
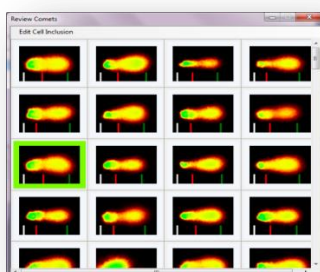
CometChip® System

CometChip® System is a high-throughput platform, shown to have a 10,000% increase in throughput compared to the original comet assay. It allows for simultaneously treating and measuring DNA damage induced by different treatments, or among different cell types on a single slide. The spatially encoded micropores ensures no overlapping of comets allowing for easy and fast analysis.



Comet Analysis Software

Breakthrough software designed to improve, standardize, and make comet analysis easier. Can be used with high throughput CometChip® and original CometAssay®. Compatible with 4X images, which allows for rapid screening (up to 400 comets in seconds when analyzing CometChip®).





CometAssay® Electrophoresis System:

The electrophoresis system, manufactured with chemically resistant materials, now comes with reusable coolant that easily slides into the cooling chamber replacing the previous water chamber. The system enables users to consistently achieve highly reproducible results for both alkaline and neutral comet assays by solving problematic variables associated with temperature, distance between electrodes and buffer height.

Features:

- Maintains constant buffer temperature with reusable cooling pack.
- Maintains optimal buffer level for consistent results with overlay.
- Specially designed trays accommodate 2, 3, 20 and 96 well CometSlides™ and CometChips® to maintain correct position during electrophoresis.
- Optimized for use with both CometAssay® and CometChip® Systems.
- Available with area specific power supply.

Ordering Information:

CometAssay®:	<u>4250-050-K*</u>	<u>4250-050-ESK*</u>	<u>4253-050-ESK-PS1*</u>
CometSlides™	x	x	x
LMAgarose	x	x	x
Lysis Solution	x	x	x
EDTA Buffer	x	x	x
Control Cells		x	x
Electrophoresis Unit		x	x
Power Supply			x

*Available with 2, 20, or 96 well slide kits

CometChip®:	4250-096-K	4260-096-CSK	4260-096-ESK	4260-096-ESK-PSI
CometChip®, 30 micron	x	x	x	x
LMAgarose	x	x	x	x
Lysis Solution	x	x	x	x
96-Well CometChip® System		x	x	x
Electrophoresis Unit			x	x
Power Supply				x

INQUIRE ABOUT BUNDLE DISCOUNTS!

Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Phone: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800 • Fax: 301-560-4973 • Email: info@trevigen.com • http://www.trevigen.com