



New, Innovative Comet Assay Electrophoresis Design

As the leading provider in standardized Comet Assay system and reagents, Trevigen announces the release of the second generation CometAssay® Electrophoresis System (ES) (4250-050-ES), designed for enhanced durability and ease of use while retaining the same electrophoresis run parameters as the original unit. 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 coolant.
- Maintains optimal buffer level for consistent results with overlay.
- Specially designed trays accommodate 2, 3, 20 and 96 well slides and maintain correct position during electrophoresis.
- Optimized for use with CometAssay Kits and CometAssay Control Cells.
- Available with area specific power supply.

Ordering Information:

Name	Catalog Number	Size	Price
CometAssay® Electrophoresis System II	4250-050-ES	1 unit	\$1,995

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>