

# TREVIGEN® Product Data

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

## CometAssay™ HT Slide

**Catalog #:** 4252-040-01  
4252-200-01  
4252-500-01  
4252-02K-01

**Size:** 2 slides  
10 slides  
25 slides  
100 slides

**Description:** 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. As a result of the migration, comets are formed and the amount of DNA in the comet tail as well as the distance of migration can be a measure of damage. Trevigen's CometSlide™ HT (High Throughput) provides a convenient higher throughput, simple and reliable sample preparation, processing, analysis and storage format for the comet assay.

**Physical State:** Sample-ready pretreated glass slides. 3 x 2 inches (75 mm X 50 mm)

**Storage:** Store at room temperature with desiccant.

**Applications:** For use with Trevigen's CometAssay™ and FLARE assay reagents and protocol.

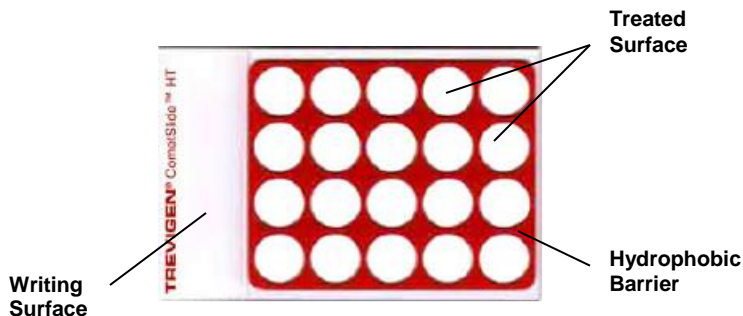


Figure 1. CometSlide™ HT features

*CometSlide™ HT*

Catalog #:

Store at Room Temperature

**TREVIGEN®**

1-800-873-8443