

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

CometAssay™ HT Slide

Catalog #: 4252-020-01	Size: 1 slide
4252-040-01	2 slides
4252-200-01	10 slides
4252-400-01	20 slides
4252-02K-01	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.

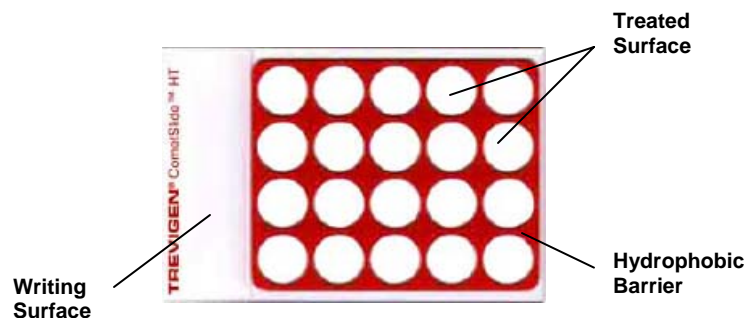


Figure 1. CometSlide™ HT features

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

CometAssay™ HT Slide

Catalog #: 4252-020-01	Size: 1 slide
4252-040-01	2 slides
4252-200-01	10 slides
4252-400-01	20 slides
4252-02K-01	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.

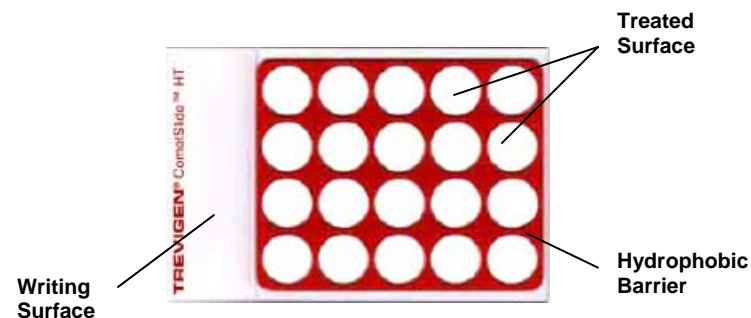


Figure 1. CometSlide™ HT features

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

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™ HT
Catalog #: 4252-02K-01
Store at Room Temperature
TREVIGEN®
1-800-873-8443

CometSlide™ HT
Catalog #: 4252-02K-01
Store at Room Temperature
TREVIGEN®
1-800-873-8443