



# TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

## Neutral CometAssay® Control Cells

### For the standardization of comet assays run under neutral conditions

Trevigen's Neutral CometAssay® Control Cells are a set of cell preparations containing different levels of DNA damage to be used as controls with Trevigen's CometAssay® Kits. When performing the neutral Comet Assay®, the four control cell populations show incremental increases in tail moment. The healthy cell population (NC0) was treated with Bleomycin under various conditions to increase the amount of damage in populations NC1, NC2 and NC3, respectively. Bleomycin complexed with metal ions induces double-stranded and single-stranded DNA breaks by a mechanism involving free radicals. These cryopreserved cells are designed to act as controls to standardize and compare comet assay results under neutral electrophoresis conditions between individual users and laboratories.

#### Ordering Information:

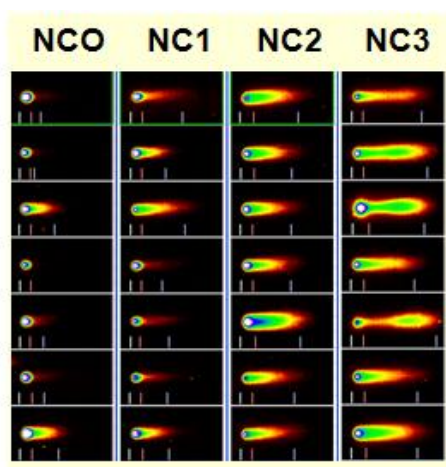
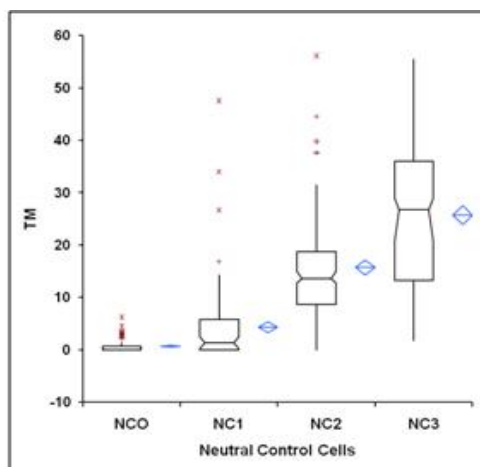
Description	Size	Catalog #
Neutral CometAssay® Control Cells	1 Set (10 Assays)	4257-010-NC

#### Kit Contents:

Description	Size	Component #
Healthy Cells	500 µl	4257-010-NC0
Treated Cells – level 1	500 µl	4257-010-NC1
Treated Cells – level 2	500 µl	4257-010-NC2
Treated Cells – level 3	500 µl	4257-010-NC3

#### Quality Control Specifications:

To ensure that our control cells allow for standardization of your comet assay results, a one-way Analysis of Variance (ANOVA) is performed to test for a significant difference between means of the treated control cells (NC1, NC2 and NC3).



Sample Images



# TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

## Neutral CometAssay® Control Cells

### Features & Benefits:

Features	Benefits
Only available set of qualified control cells for use with the neutral CometAssay®	Standardize and validate neutral electrophoresis methods between individual users and laboratories.
Includes four populations: one population of healthy cells and three populations with increasing amounts of DNA damage	Provides a range of DNA damage for comparison. Damaged populations show statistically distinct differences.
Cells are cryopreserved	After initial thaw, aliquot and refreeze, the cells remain stable in liquid nitrogen for long term storage.
Qualified for neutral gel electrophoresis	Standardized with Trevigen's CometAssay® Electrophoresis System (cat# 4250-050-ES) designed to eliminate known causes of assay variability.

### CometAssay® Product Portfolio:

Description	Size	Catalog #
CometAssay® Electrophoresis System	1 System	4250-050-ES
CometAssay® Electrophoresis Starter Kit (includes CometAssay® Kit, Control Cells (alkaline) & ES)	Starter Kit	4250-050-ESK
CometAssay® Kit	1 Kit	4250-050-K
CometAssay® Silver Kit	1 Kit	4251-050-K
CometAssay® Silver Staining Kit	1 Kit	4254-200-K
CometAssay® HT Sample Kit	1 Kit	4252-040-K
CometAssay® Kit 96 Wells	1 Kit	4253-096-K
CometSlide™ (2-well slide)	25 Slides	4250-050-03
CometSlide™ (2-well slide)	100 Slides	4250-200-03
CometAssay® HT Slide (20 well slide)	25 Slides	4252-500-01
CometAssay® HT Slide (20 well slide)	100 Slides	4252-02K-01
96-Well CometSlide™ (96 well slide)	10 Slides	4253-960-03
96-Well CometSlide™ (96 well slide)	20 Slides	4253-02K-03
96-Well CometSlide™ (96 well slide)	100 Slides	4253-10K-03
CometAssay® Control Cells (alkaline)	1 Set (10 Assays)	4256-010-NC
Neutral CometAssay® Control Cells	1 Set (10 Assays)	4257-010-NC

©2011 CometSlide is a trademark of Trevigen, Inc. Trevigen and CometAssay are registered trademarks of Trevigen, Inc.

Trevigen, Inc.  
8405 Helgerman Court  
Gaithersburg, Maryland 20877  
1-800-873-8443