

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Neutral CometAssay® Control Cells

Catalog #: 4257-010-NC

Lot#: 38986C17

Components

4257-010-NC0
 4257-010-NC1
 4257-010-NC2
 4257-010-NC3

Component Lot

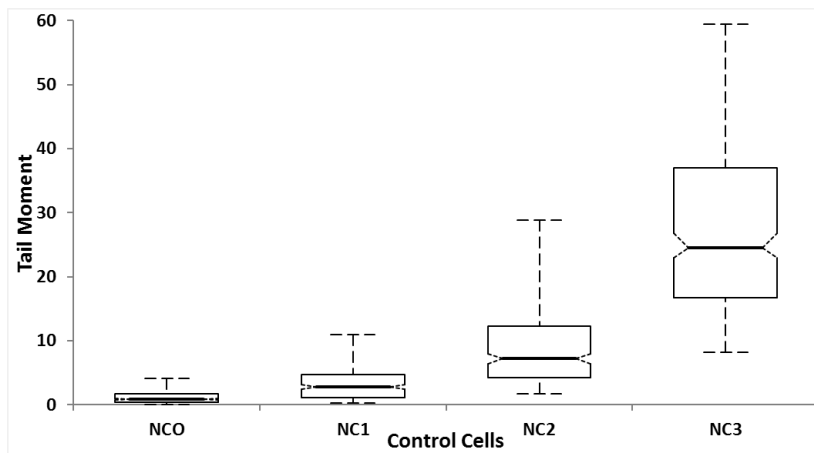
38982C17
 38983C17
 38984C17
 38985C17

Description: Trevigen's Neutral CometAssay® Control Cells are a set of asynchronous cell populations comprised of a healthy population (NC0) and cells treated with Bleomycin to increase the amount of DNA damage in populations NC1, NC2, and NC3, respectively.

Storage Conditions: Neutral CometAssay® Control Cells are stored using a Liquid Nitrogen Storage System.

Assay Conditions: These cell preparations containing different levels of DNA damage are to be used as controls for comet assays run under neutral conditions with Trevigen's CometAssay® Kits. Lot specific data is generated to calculate Tail Moment for each Control Cell population.

Lot Specific Data: Box-Whisker plot of Control Cells: Tail Moment



TM by Control Cells	N	Mean	80% CI	SE	SD	Median	80% CI	IQR
NCO	153	1.393	0.359 to 2.428	0.8064	2.066	0.850	0.730 to 0.980	1.343
NC1	206	4.001	3.110 to 4.893	0.6950	5.374	2.795	2.390 to 3.210	3.636
NC2	191	10.391	9.466 to 11.317	0.7218	11.695	7.230	6.420 to 8.020	7.993
NC3	169	28.321	27.337 to 29.305	0.7673	15.155	24.510	22.970 to 26.800	20.287

©2017 Trevigen, Inc. All rights reserved. Trevigen, and CometAssay are registered trademarks, and CometSlide is a trademark of Trevigen, Inc. E3/08/17v1

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@trevigen.com • www.trevigen.com

Neutral CometAssay[®] Control Cells
Catalog #: 4257-010-NC
Store in Liquid Nitrogen
TREVIGEN[®]
1-800-873-8443