

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Alkaline CometAssay® Control Cells **Catalog #:** 4256-010-CC **Lot #:** 38594A17

Components

4256-010-CC0
 4256-010-CC1
 4256-010-CC2
 4256-010-CC3

Component Lot #

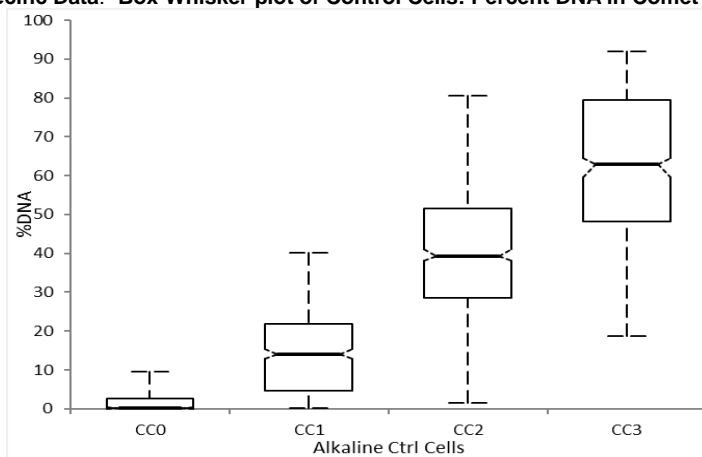
38583A17
 38584A17
 38585A17
 38586A17

Description: Trevigen's Alkaline CometAssay® Control Cells are asynchronous cell populations composed of a healthy population (CC0) and cells treated with Etoposide to increase the amount of damage in populations CC1, CC2, and CC3, respectively.

Storage Conditions: 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 with Trevigen's CometAssay® Kits. Lot specific data is generated to calculate percent DNA in tail for each Control Cell population.

Lot Specific Data: Box-Whisker plot of Control Cells: Percent DNA in Comet Tail



% DNA in Tail by Alkaline Ctrl Cells	N	Mean	Mean SE	SD	80% CI*	Median	IQR	80% CI
CC0	490	2.420	0.7642	5.948	1.440 to 3.400	0.250	2.560	0.240 to 0.290
CC1	334	15.610	0.9256	13.573	14.423 to 16.796	14.020	17.233	12.810 to 15.280
CC2	462	40.155	0.7870	20.703	39.146 to 41.164	39.260	23.145	38.150 to 40.970
CC3	391	61.231	0.8555	22.787	60.134 to 62.328	62.870	31.308	59.680 to 64.500

©2017 Trevigen, Inc. All rights reserved. Trevigen and CometAssay are registered trademarks, and CometSlide is a trademark of Trevigen, Inc.

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

CometAssay® Control Cells
Catalog #: 4256-010-CC
Lot #: 38594A17
Store in Liquid Nitrogen
TREVIGEN®
1-800-873-8443