

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Alkaline CometAssay® Control Cells Catalog #: 4256-010-CC Lot #: 40080G17

Components

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

Component Lot #

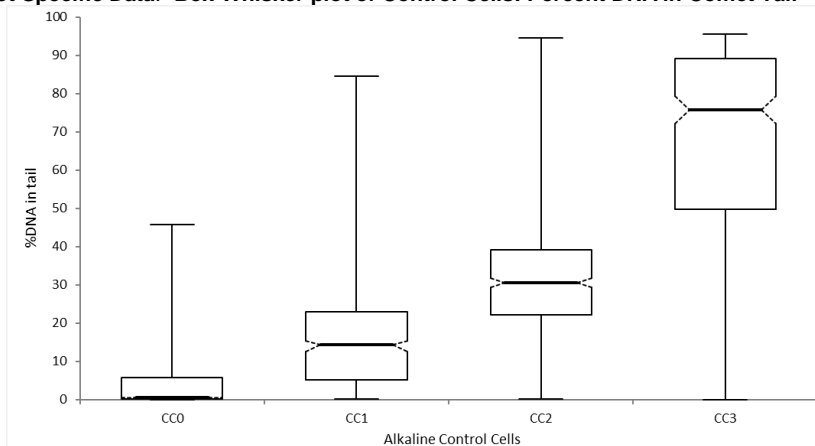
40065G17
 40066G17
 40067G17
 40068G17

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 Control Cells	N	Mean	Mean SE*	SD	80% CI†	Median	IQR	80% CI
CC0	237	3.954	1.0583	6.473	2.597 to 5.311	0.440	5.590	0.370 to 0.510
CC1	250	15.393	1.0304	12.770	14.071 to 16.714	14.240	17.767	12.770 to 15.250
CC2	238	31.966	1.0561	15.102	30.612 to 33.320	30.625	17.018	29.340 to 31.660
CC3	198	67.145	1.1578	26.601	65.660 to 68.630	75.655	39.419	72.240 to 79.280

©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 #: 40080G17
Store in Liquid Nitrogen
TREVIGEN®
1-800-873-8443