

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

CometAssay® Alkaline Control Cells Catalog #: 4256-010-CC Lot #: 37035G16

Components

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

Component Lot #

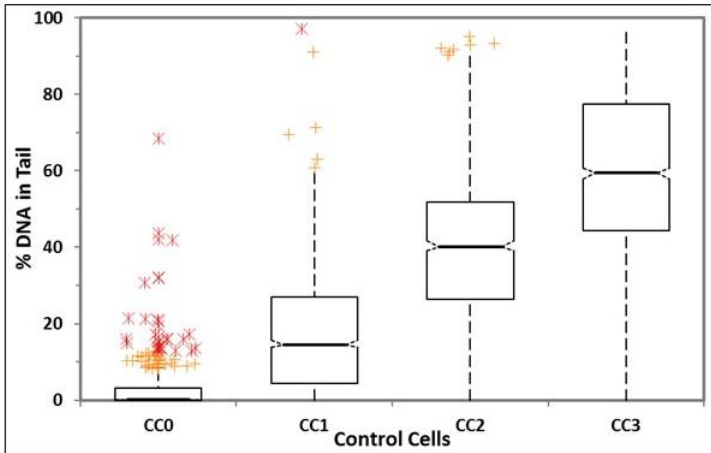
32553B15
 32554B15
 37018G16
 32556B15

Description: Trevigen’s CometAssay® Alkaline 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 Control Cells	N	Mean	SE	SD	80% CI	Median	IQR	80% CI
CC0	537	2.826	0.7213	9.186	1.902 to 3.751	0.150	3.230	0.130 to 0.170
CC1	669	17.869	0.6462	17.611	17.040 to 18.697	14.550	22.493	13.810 to 15.590
CC2	599	39.353	0.6829	19.608	38.478 to 40.229	40.180	25.552	39.030 to 41.650
CC3	430	58.993	0.8060	18.883	57.959 to 60.026	59.405	33.093	57.890 to 60.760

©2016 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 #: 37035G16
Store in Liquid Nitrogen
TREVIGEN®
1-800-873-8443