

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Alkaline CometAssay® Control Cells Catalog #: 4256-010-CC Lot #: 41002L17

Components

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

Component Lot #

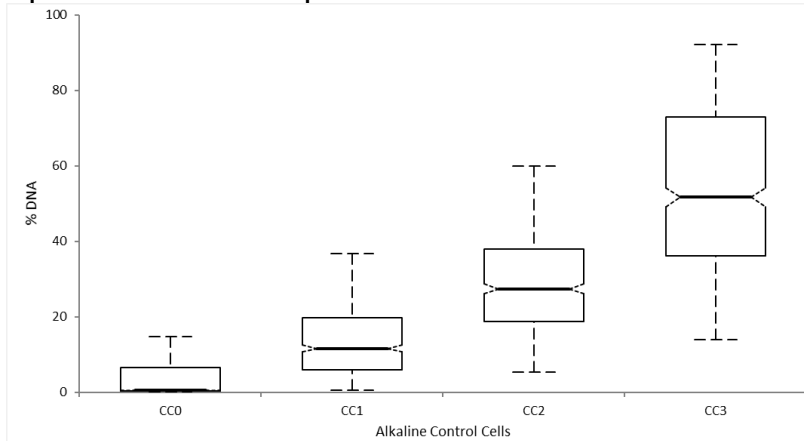
40888L17
 40889L17
 40890L17
 40891L17

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 by Alkaline Control Cells	N	Mean	Mean SE	SD	80% CI	Median	IQR
CC0	378	3.735	0.2785	5.414	0.420 to 0.550	0.440	6.278
CC1	377	14.105	0.6156	11.953	10.650 to 12.530	11.590	13.843
CC2	372	29.343	0.8538	16.467	26.160 to 28.810	27.385	19.114
CC3	455	53.573	1.1254	24.006	49.070 to 54.180	51.700	36.967

©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 #: 41002L17
Store in Liquid Nitrogen
TREVIGEN®
1-800-873-8443