

# Trevigen Cell Assays

*Trevigen Cell Assays, (TCA) a division of Trevigen, Inc, was established in 2008 to conduct contract research work for the pharmaceutical, biotechnology, government and academic segments of the research market. TCA specializes in designing and conducting assays for lead compound and genotoxic screening based on DNA damage and repair and cancer cell behavior.*

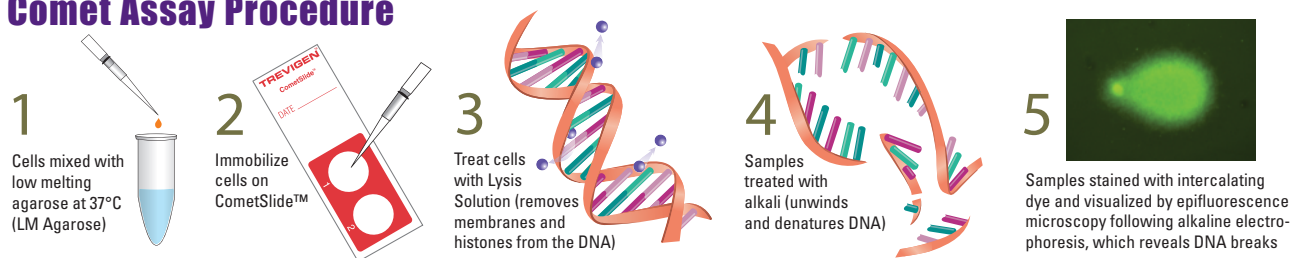
**TREVIGEN®**

COMET ASSAY

## Comet Assay Service

Trevigen is the sole provider of a standardized CometAssay® system for the direct detection of DNA Damage. The CometAssay® is a single cell gel electrophoresis assay which can be used to quantitate DNA Damage from in vivo or in vitro samples. The CometAssay® measures double strand breaks (DSBs), single strand breaks (SSBs), alkali labile sites, oxidative DNA base damage, DNA-DNA/DNA-protein/DNA-Drug crosslinking and DNA repair. At your request, the scientists at TCA will collaborate with you to design screening studies employing the CometAssay® tailored for your specific needs.

## Comet Assay Procedure



## Advantages

- TCA uses a unique platform of standardized comet assay kits, slides, control cells, CometAssay® ES electrophoresis system (patent applied for) and state of the art image analysis to assure consistency of results with successive experiments or screening studies.
- The technical team at TCA developed the reagents, slides, control cells and electrophoresis system, and is unmatched in familiarity and skill with the comet assay.
- TCA has already successfully executed multiple comet assay contracts and thus has a proven track record of high performance.
- Confidentiality is assured. TCA is ready to execute Non-Disclosure Agreements as required.

## Also Available

TCA has expanded its services to offer study design and compound screening services for:

- PARP *in vivo* Pharmacodynamic Assay II (FORM TCA 2)
- PARP/PARG Assays (FORM TCA 3)
- Cell Invasion/Migration (FORM TCA 4)

# Trevigen Cell Assays

**COMET ASSAY**

Getting started is easy. Fax the completed form to 301-560-4973; or email us at [TCA@trevigen.com](mailto:TCA@trevigen.com); or complete the form online at [www.trevigen.com](http://www.trevigen.com) with the information that we need in order to set up your screening service.

## Quotation Request Form - Comet Assay

First and Last Name \_\_\_\_\_

Email Address \_\_\_\_\_

Company \_\_\_\_\_

Street Address \_\_\_\_\_

City \_\_\_\_\_ State/Province \_\_\_\_\_

Zip/Postal Code \_\_\_\_\_ Country \_\_\_\_\_

Telephone Number \_\_\_\_\_ Fax Number \_\_\_\_\_

How many samples do you have to screen? \_\_\_\_\_

How many replicates are required? \_\_\_\_\_

Are you providing the samples or will treatment be done by Trevigen Cell Assays? \_\_\_\_\_

A) How many compounds and concentrations do you have to screen? \_\_\_\_\_

B) How many time points? \_\_\_\_\_ What is the treatment duration? \_\_\_\_\_

C) What types of cells are required? \_\_\_\_\_

D) Will you be providing the cells? \_\_\_\_\_

What is the desired reporting format? \_\_\_\_\_

How soon is the data required? \_\_\_\_\_

Are there any other screening parameters or special conditions that you require? \_\_\_\_\_

### Compound Handling Instructions

What compound(s) are you screening? \_\_\_\_\_

Will you be providing the compound(s)? \_\_\_\_\_

If you are not providing the compound(s) where can they be purchased? \_\_\_\_\_

Is the compound toxic? If yes, are MSDS available? \_\_\_\_\_

What storage conditions are required? \_\_\_\_\_

What diluent is required? \_\_\_\_\_

Upon receipt, a TCA senior scientist will contact you to go over the desired work and discuss options as appropriate. A proposal and cost will then be prepared. The proposal will include the turn around time and the agreed upon reporting format.