



Trevigen Cell Assay Service

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 medical research market. TCA specializes in designing and conducting assays for lead compounds and genotoxic screening based on DNA damage and repair as well as cancer cell behavior.

PROCESS:

- A Project Manager and a Senior Scientist is assigned for every project
- A detailed Protocol and Schedule is prepared before the work begins
- Site training for sample collection is available if applicable
- Observations, comments, and updates are sent by e-mail during the study
- A Final Report with raw data and basic Statistical/Graphical analysis is provided.

ADVANTAGES:

- The technical team at TCA developed the assays and is unmatched in skill and knowledge
- Confidentiality is assured. A Non-Disclosure Agreement can be executed if requested

TCA Service:	Assay:	Application:
CometAssay® Cat# 4250-050-TCA	Rapid analysis of DNA fragmentation associated with DNA damage.	<ul style="list-style-type: none"> • Detect and quantitate DNA damage • Follow DNA repair
PARP Assay Cat# 4676-096-TCA Cat# 4677-096-TCA	Measurement of biotinylated poly(ADP-ribose) incorporation onto histone proteins.	<ul style="list-style-type: none"> • Assay inhibitors and activators of PARP activity • Determination of IC₅₀ values for PARP inhibitors
PARG Assay Cat# 4682-096-TCA Cat# 4683-096-TCA	Measurement of the loss of biotinylated PAR from histones.	<ul style="list-style-type: none"> • Identify inhibitors and activators of PARG activity • Determination of IC₅₀ values for PARG inhibitors
Pharmacodynamic PARP Assay Cat# 4520-096-TCA	Measurement of new PAR levels in cellular extracts and tumor lysates.	<ul style="list-style-type: none"> • Quantification of PAR in peripheral blood mononuclear cells, tissue culture cells, and tumor lysates from different tissues, organs and xenografts • Monitoring the efficacy of PARP inhibitors on PAR formation <i>in vivo</i>
Cell Invasion Assay Cat# 3455-024-TCA Cat# 3465-096-TCA	Analysis of responses of chemokines, toxins, drugs and other analytes of interest.	Screening for compounds that influence cellular invasion through <ul style="list-style-type: none"> • Extracellular matrices, • Angiogenesis • Immune response • Tumor cell metastasis
Tube Formation Assay Cat# 3470-096-TCA	Investigation of prospective angiogenesis inhibitors.	<ul style="list-style-type: none"> • Detection of inducers and inhibitors of endothelial cell tube formation

*Please call 1-800-873-8443 or visit us at www.trevigen.com and click on the services tab for more information and custom quotation forms.

Trevigen Cell Assay Service