

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.

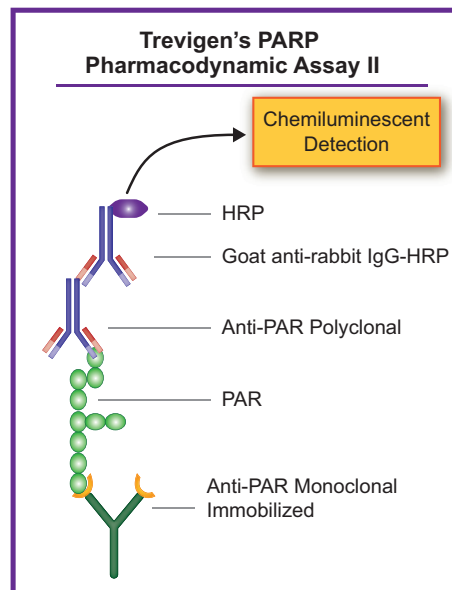
TREVIGEN®

PARP PHARMACODYNAMIC ASSAY

HT PARP *in vivo* Pharmacodynamic Assay

The PDA II kit is for RESEARCH USE ONLY

PARP catalyzes the NAD⁺-dependent addition of poly (ADP-ribose) (PAR) onto itself and adjacent nuclear proteins. This enzyme is a therapeutic target for BRCA1 and BRCA2 associated breast cancers. To address the need to monitor PARP activity among different individuals and within cells. Trevigen's improved and validated HT PARP *in vivo* Pharmacodynamic Assay II measures net PAR levels in tissue or cellular extracts and has been used to document differences in PAR levels among tumor lysates, organs and xenografts. The HT PARP *in vivo* Pharmacodynamic Assay II employs a 96 well plate, pre-coated with Trevigen's monoclonal PAR antibody as the capture agent, and anti-PAR polyclonal rabbit antibody as the detecting agent.



Assay Design

Step 1: Immobilized PAR mAb captures cellular PAR and PAR attached to proteins in prepared lysates.

Step 2: Binding of PAR polyclonal detecting Ab to capture PAR.

Step 3: Measure captured PAR via binding of goat anti-rabbit IgG-HR with chemiluminescent detection.

• Light output (Signal) correlates with the amount of cellular PAR.

	Catalog No.	Application(s)	Detection Method	Test Sample	Sensitivity
PARP <i>in vivo</i> Pharmacodynamic Assay II • Chemiluminescence Detection	4520-096-K	PARP Inhibition in vivo PAR Levels in Lysates	PAR Ab In vivo PARylation	PMBC Lysates Tissue Lysates	2-1000 pg/ml PAR

Trevigen Cell Assays

FORM TCA 2 DATE

PARP PHARMACODYNAMIC ASSAY

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

Quotation Request Form - PDAll Assay

First and Last Name _____
Email Address _____
Company _____
Street Address _____
City _____ State/Province _____
Zip/Postal Code _____ Country _____
Telephone Number _____ Fax Number _____

General Questions

What is the desired reporting format? _____ How soon is the data required? _____
How many samples will be tested? _____ How many replicates are required? _____
Are there any other screening parameters or special conditions that you require? _____
What type of samples are being tested? _____

If Clinical Research Samples:

Are the samples part of a clinical trial? _____
If yes, will coordination with different clinical sites be required? _____
Do you require training in sample preparation? _____

If Drug Discovery Samples:

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

If Treatment Is Being Performed By Trevigen Cell Assays

How many compounds and concentration range to be tested do you have to screen? _____
How many time points? _____ What is the treatment duration? _____
What type of cells will be used? _____
Will you be providing the cells? _____
Do the cells require any special handling? _____
What type of damaging agent is required? _____

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 is solubility of compound? _____

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.

Trevigen, Inc.
8405 Helgerman Court
Gaithersburg, Maryland 20877
info@trevigen.com
1-800-873-8443 • www.trevigen.com