

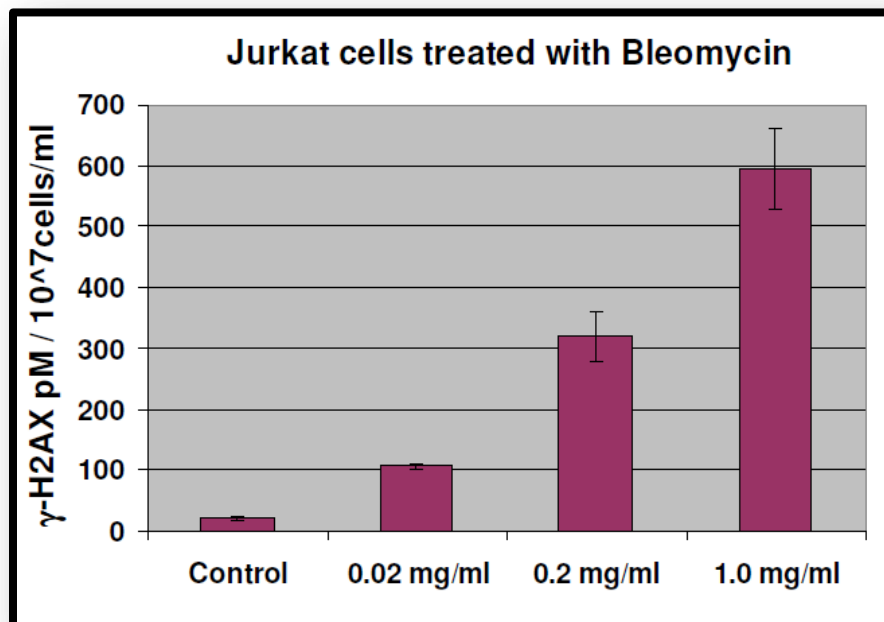


Trevigen®

Better Products. Better Results.

γ -H2AX Pharmacodynamic ELISA Kit

Trevigen's γ -H2AX Pharmacodynamic Assay is the first commercially available kit for the study of double strand DNA breaks through the detection of γ -H2AX, a phosphorylated histone historically proven as a highly specific and sensitive molecular marker for double strand DNA break detection. The assay documents γ -H2AX levels in peripheral blood mononuclear cells, cultured cells and tissue biopsies, and is available as a complete reagent kit with chemiluminescent detection.



γ -H2AX values from Jurkat Cells treated with Bleomycin. Jurkat cells treated at 5×10^6 cells/ml with 0.02 mg/ml, 0.2 mg/ml and 1.0 mg/ml of Bleomycin respectively.

FEATURES:

- Chemiluminescent, non-radioactive ELISA
- Pre-coated 96 well capture antibody plate
- Dynamic range from 10 pM to 800pM
- Sensitivity with 5 pM of γ -H2AX

APPLICATIONS: quantitates γ -H2AX levels in PBMC, cultured cells and tissue biopsies.

KIT COMPONENTS:

96-stripwell pre-coated plate with 3 film sealers
 γ -H2AX Standard
Assay Buffer
H2AX IgM Detecting Antibody
Goat anti-Mouse IgM HRP Conjugate

Cell Lysis Reagent
Jurkat Cell Lysate Control
25X Wash Buffer
PeroxyGlow™ A
PeroxyGlow™ B

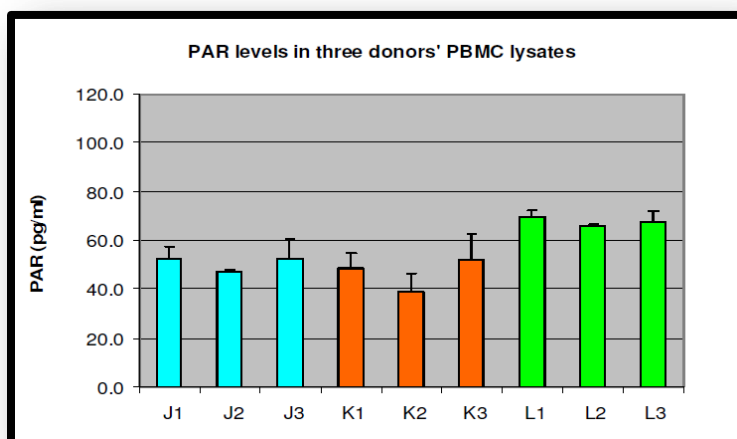
ORDERING INFORMATION:

Name	Catalog Number	Size	Price
γ -H2AX Pharmacodynamic Assay	4418-096-K	96 samples	\$1,250



PARP in vivo Pharmacodynamic ELISA Kit II

Trevigen's improved and validated 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 assay 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.



Baseline PAR levels in Peripheral Blood Mononucleocytes (PBMCs) from normal donors (J, K, and L) expressed in terms of pg/ml per 1×10^7 cells/ml. PBMC lysates were made on three days and each was assayed in triplicates. The means and standard deviations of each determination are shown.

FEATURES:

- Validated ELISA assay that measures drug action on PARP in both in vivo and in vitro setting
- Pre-coated 96 well capture antibody plates
- High signal to noise ratio
- Detection sensitivity of 2 pg/ml of PAR
- Broad linear dynamic range to 1,000 pg/ml
- Reduced inter-assay variability

APPLICATIONS:

- Quantitation 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 in vivo.
- Verifying observations of enhanced cancer cell cytotoxicity arising from PARP inhibitor/anticancer drug combination therapy.
- Facilitating development of PARP and PARG targeted therapeutics.

KIT COMPONENTS:

PAR Standard and Sample Buffer
PAR Polyclonal Detecting Ab and Diluent
Goat anti-Rabbit IgG-HRP
Detection Reagents
Cell Lysis Reagent and 20% SDS

DNase I and Mg Cation
Jurkat Cell Lysate Standard Controls
(Low, Medium, High)
Pre-coated 96-stripwell plate and seal

ORDERING INFORMATION:

Name	Catalog Number	Size	Price
PARP in vivo Pharmacodynamic Assay II	4520-096-K	96 samples	\$1,250

Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Phone: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800 • Fax: 301-560-4973 • Email: info@trevigen.com • <http://www.trevigen.com>