

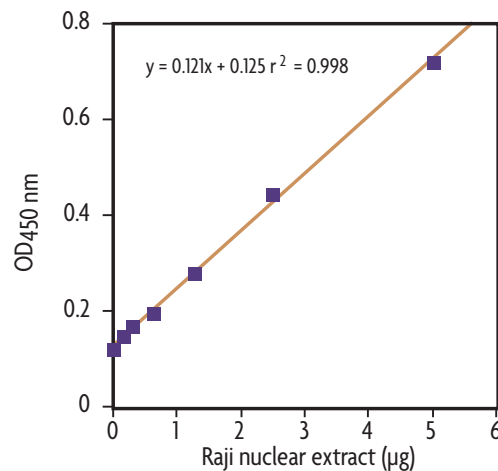
DNA Repair Protein ELISAs

rapid, sensitive analysis of GTBP, Ku and RPA protein activity

DNA repair proteins maintain genome integrity by recognizing, binding to, and repairing damaged DNA. Deficiencies in their activity have been linked to many different pathological diseases, including cancer. To facilitate the study of DNA repair protein activity in tissue and cell culture extracts, Active Motif developed its DNA Repair Protein Kits. These kits are ELISA-based assays that include a 96-well plate in which multiple copies of a specific oligonucleotide that contains DNA damage have been immobilized. When extract is added, the repair protein of interest binds to the oligonucleotide. Incubation with primary antibody that is specific for the repair protein being studied, and with secondary HRP-conjugated antibody and developing solution provides an easily quantified colorimetric readout.

The DNA Repair Protein ELISA advantage

- Sensitive, colorimetric readout easily quantified by spectrophotometry
- ELISA format eliminates gels, blotting and radioactivity
- From cell extract to completed assay in less than 5 hours
- Up to 10-fold greater sensitivity than gelshift assays
- Ability to assay both cells and tissues
- High-throughput compatible



Monitoring Ku70 DNA binding activity using the Ku70/86 DNA Repair Kit.

Nuclear extract from untreated Raji cells is tested for Ku activity using the Ku70/86 DNA Repair Protein Kit.

CONTENTS

One or five 96-well plates with plate sealer(s), Primary Antibody(ies), HRP-conjugated secondary antibody, positive control cell extract, competitor oligos, Lysis, Binding and Wash Buffers, Protease Inhibitor, DTT and Developing Solutions.

Product	Format	Catalog No.
GTBP DNA Repair Protein Kit	1 x 96 rxns	51096
	5 x 96 rxns	51596
Ku70/86 DNA Repair Protein Kit	1 x 96 rxns	51196
	5 x 96 rxns	51696
RPA DNA Repair Protein Kit	1 x 96 rxns	51296
	5 x 96 rxns	51796