

Nuclear Receptors

kits and reagents for studying nuclear receptor protein activity

Nuclear Receptors (NRs) are important regulators of genomic activity; inappropriate signaling by NRs is associated with many different disease states, including cancer, asthma, infertility, osteoporosis and arthritis. As such, NRs have proven to be promising drug targets. With Active Motif's complete line of kits and reagents for studying ER, GR, PPAR, VDR and more, you now have access to a variety of products that will help you effectively study NRs. We offer TransAM™ ELISAs for monitoring the DNA binding activity of NRs, NR Peptide ELISAs for monitoring ligand-activated NRs, Sandwich ELISAs to examine total amounts of NR proteins, as well as a vast line of NR antibodies and recombinant proteins suitable for a variety of applications such as ChIP, Western blot and EMSA.

The Nuclear Receptor product line

- Complete solution for studying nuclear receptors
- Monitor DNA binding activity of NRs with TransAM™ ELISAs (Figure 1)
- Compare agonist/antagonist effects using NR Peptide ELISAs (Figure 2)
- Monitor total NR protein levels using NR Sandwich ELISAs
- Western blot with confidence using highly characterized antibodies coupled with recombinant proteins as positive controls
- Perform traditional EMSAs using Nushift™ and Gelshift™ Kits

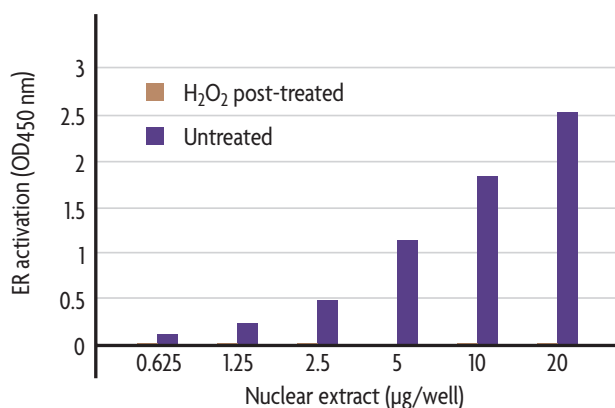


Figure 1: Monitoring ERα DNA binding activity using TransAM ER.

Different amounts of untreated and H₂O₂ post-treated nuclear extracts from MCF-7 cells are tested for ER activity using the TransAM ER Kit.

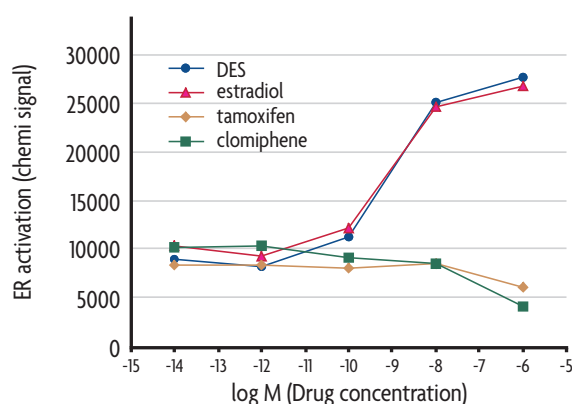


Figure 2: ERα agonism/antagonism dose-response curve of test compounds.

MCF-7 nuclear extracts are incubated in the presence of DES, estradiol, tamoxifen and clomiphene in wells of the NR Peptide Binding ELISA. Only ligand-activated ER binds to the Capture Peptide on the plate.

TransAM™ DNA Binding ELISAs

Product	Format	Catalog No.
TransAM™ ER	1 x 96 rxns	41396
	5 x 96 rxns	41996
TransAM™ GR	1 x 96 rxns	45496
	5 x 96 rxns	45996
TransAM™ PPAR γ	1 x 96 rxns	40196
	5 x 96 rxns	40696

Peptide ELISAs

Product	Format	Catalog No.
NR Peptide ER α	1 x 96 rxns	49096
	5 x 96 rxns	49596
NR Peptide ER α Chemi	1 x 96 rxns	49097
	5 x 96 rxns	49597

Sandwich ELISAs

Product	Format	Catalog No.
NR Sandwich AR	1 x 96 rxns	49196
	5 x 96 rxns	49696
NR Sandwich ER α	1 x 96 rxns	49296
	5 x 96 rxns	49796
NR Sandwich Total PR	1 x 96 rxns	49396
	5 x 96 rxns	49896
NR Sandwich PR B	1 x 96 rxns	49496
	5 x 96 rxns	49996

Nushift™ and Gelshift™ Kits

Product	Format	Catalog No.
Nushift™ ER α	17 rxns	37050
Nushift™ PPAR γ	17 rxns	37034
Nushift™ VDR	17 rxns	37042
Gelshift™ ER	20 rxns	37309
Gelshift™ HNF-4	20 rxns	37312
Gelshift™ PPAR	20 rxns	37321

Recombinant Proteins

Product	Format	Catalog No.
Recombinant ER Protein	4,000 ng	31119
Recombinant FXR Protein	10,000 ng	31120
Recombinant GR Protein	5,000 ng	31121
Recombinant LXR α Protein	10,000 ng	31122
Recombinant LXR β Protein	10,000 ng	31123
Recombinant PPAR α Protein	10,000 ng	31125
Recombinant PPAR β (δ) Protein	10,000 ng	31126
Recombinant PPAR γ Protein	10,000 ng	31127
Recombinant RAR- α Protein	5,000 ng	31130
Recombinant RAR- β Protein	5,000 ng	31131
Recombinant RAR- γ Protein	5,000 ng	31132
Recombinant RXR- α Protein	10,000 ng	31133
Recombinant RXR- β Protein	5,000 ng	31134
Recombinant RXR-LBD Protein	10,000 ng	31135

Antibodies

Product	Format	Catalog No.
ER α mAb for WB	100 μ l	39510
ER α pAb for WB	100 μ l	39021
ER α pAb for EMSA	17 rxns	39316
GR pAb for WB	100 μ l	39027
HNF-4 pAb for WB	100 μ g	39076
PPAR α pAb for WB	100 μ l	39045
PPAR γ pAb for EMSA & CHIP	17 rxns	39338
PXR pAb for WB	100 μ l	39048
RAR- β pAb for WB	100 μ l	39049
RAR- β 2 pAb for WB	100 μ l	39050
RAR- γ pAb for WB	100 μ l	39051
RXR- α pAb for WB	100 μ l	39054
RXR- β pAb for WB	100 μ l	39055
VDR pAb for WB	100 μ l	39069
VDR pAb for EMSA	17 rxns	39344