

# Nuclear Extract Kit

high yields of specifically segregated extracts

Active Motif's Nuclear Extract Kit is ideal for the preparation of nuclear, cytoplasmic and whole-cell extracts from mammalian cells and tissues. The resultant high-quality extracts may be used with TransAM™ Kits or in gelshift assays, Western blots, DNA footprinting or as a starting point for transcription factor purification.

The Nuclear Extract Kit eliminates the need to optimize reagents and ensures consistently high yields. The detailed protocol helps ensure that your extract is not contaminated with proteins from other cellular compartments (Figure 1). A standardized extract procedure also eliminates variability from your experiments, so you'll get reproducible results.

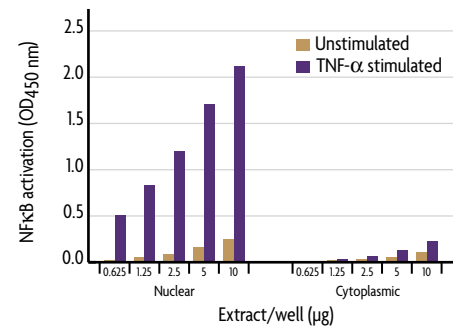
Product	Format	Catalog No.
Nuclear Extract Kit	100 rxns	40010
	400 rxns	40410

## CONTENTS

Lysis Buffer, Dithiothreitol, Protease Inhibitor Cocktail, 10X PBS, Phosphatase Inhibitors, 10X Hypotonic Buffer and Detergent. Sufficient reagents are included for 100 or 400 extractions from a 100 mm dish.

## NUCLEAR EXTRACT KIT ADVANTAGES

- Complete kit contains all required reagents
- No need to optimize your own procedure
- Quality-controlled reagents ensure reproducibility



**Figure 1: Specific extraction of nuclear and cytoplasmic extracts.**

Nuclear and cytoplasmic extracts were prepared using the Nuclear Extract Kit from HeLa samples unstimulated or stimulated with TNF- $\alpha$  for 30 minutes and assayed using the TransAM™ NF $\kappa$ B p50 Kit. Because activated NF $\kappa$ B is translocated to the nucleus, only nuclear extract from stimulated cells should contain activated NF $\kappa$ B.