

# KinaseSTAR™ JNK Activity Screening Kit

(Catalog #K430-40; 40 assays; Store kit at -20°C)

## I. Introduction:

c-Jun N-terminal kinase (JNK) is one of the several main MAP kinase groups identified in mammals. Recent evidences suggest that activation of JNK may play an important role in neuronal apoptosis and other physiological and pathological processes. The JNK Activity screening kit utilizes an N-terminal c-Jun (1-79) fusion protein bound to glutathion sepharose beads to selectively “pull down” JNK from cell lysates. After wash to remove nonspecifically bound proteins, the kinase reaction is then carried out in the presence of cold ATP. c-Jun phosphorylation is then measured by Western blot analysis using a phospho-c-Jun specific antibody.

## II. Kit Contents:

Component	K430-40	Cap Color
Kinase Extraction Buffer	80 ml	NM
c-Jun (1-79) Fusion Protein Beads	800 µl	Red
Kinase Assay Buffer	25 ml	WM
ATP (10 mM)	50 µl	Yellow
Phospho-cJun Specific Antibody	50 µl	Green

## III. JNK Activity Immunoassay Protocol

### A. Preparation of Cell Lysate:

1. Activate cells by desired methods. Concurrently incubate a negative control culture without activation. To generate a positive control, cells can be treated with 1 µg/ml of Anisomycin (Cat.# 1549-10) for 1 hr, before harvested.
2. Pellet cells (2-10 millions/assay) and wash once in 1X ice-cold PBS.
3. Lyse cells in 200 µl ice-cold JNK Extraction Buffer. Incubate on ice for 5 minutes.
4. Pellet at 13,000 rpm for 10 min at 4°C. Transfer supernatant (This is the Cell Lysate) to a new tube.
5. Assay protein concentration of the Cell Lysate. The Cell Lysate can be used immediately or freeze at -80°C for future use.

### B. “Pull Down” JNK Using c-Jun Fusion Protein:

6. For each assay, add 20 µl c-Jun Fusion Protein Beads to 200 µl Cell Lysate (~50-400 µg total protein). Incubate with gentle rocking overnight at 4°C.
7. Microcentrifuge (14,000 rpm) for 30 seconds at 4°C. Remove Supernatant. Wash pellet twice with 0.5 ml of 1X Kinase Extraction Buffer and one time with 0.5 ml Kinase Assay Buffer. Keep on ice.

### C. Kinase Assay:

8. Suspend pellet in 50 µl Kinase Assay Buffer. Add 1 µl of 10 mM ATP. Incubate for 30 minutes at 30°C.
9. Add 30 µl 3X SDS-PAGE Buffer (not provided)
10. Boil the samples for 3 minutes. Microcentrifuge for 2 minutes.
11. Load the supernatant (20 µl) on 12% SDS-PAGE. Alternatively, the supernatant may be stored at -20°C for future use.

### D. Western Immunoblotting:

12. Perform Western blot analysis using the rabbit anti-Phospho-cJun (Ser 73) Specific Antibody at 1:1000 dilutions. A 35 kDa band corresponding to the phosphorylated c-Jun protein should be detected in JNK activated samples.

## RELATED PRODUCTS:

### Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set
- Apoptosis siRNA Vectors

### Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

### Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

### Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

### Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

### Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- PLTP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

### Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

### Growth Factors and Cytokines

### Monoclonal and Polyclonal Antibodies