

## Anti-phospho-ASK1 (Ser966)

Rabbit Polyclonal Antibody

### Catalog # M13-65R

Lot # J1178-2

### Cited Applications

IHC

Suggested Dilutions:

IHC: 1:50-1:100

Ideal working dilutions for each application should be empirically determined by the investigator.

### Specificity

Recognizes the ASK1 protein phosphorylated at serine 966

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser966

### Formulation

PBS (pH 7.4) 150mM NaCl, 0.02% sodium azide and 50% glycerol.

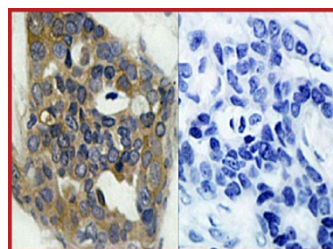
### Scientific Background

ASK1, also known as MAPKKK5, is responsible for activating MKK3, MKK4 (SEK1), and MKK6. ASK1 is activated in cells treated with tumor necrosis factor-alpha and overexpression of ASK1 leads to increased apoptosis (1). ASK1 interacts with members of the TRAF family and is activated by TRAF2 in the TNF-signaling pathway leading to the activation of MKK4 and subsequently JNK. Thus, ASK1 is a mediator of TRAF2-induced JNK activation (2).

### References

1. Ichijo, H. et al: Induction of apoptosis by ASK1, a mammalian MAPKKK that activates SAPK/JNK and p38 signaling pathways. *Science* 275: 90-94, 1997.
2. Nishitoh, H. et al: ASK1 is essential for JNK/SAPK activation by TRAF2. *Molec. Cell* 2: 389-395, 1998.

### Sample Data



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using anti-phospho-ASK1 (Ser966) antibody.

## Anti-phospho-ASK1 (Ser966)

Rabbit Polyclonal Antibody

Catalog Number

M13-65R

Specific Lot Number

J1178-2

Purification Affinity chromatography

Stability 1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**