

## Recombinant IKKε protein

**Catalog No:** 31177

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant IKKε protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.069 µg/µl.

**Background:** **IKKε – IκB Kinase-epsilon** (also called **IKKi**) is a homolog of the IKK (IκB kinase) catalytic subunits IKKα and IKKβ. IKKε is primarily expressed in immune cells and is responsible for modulation of IRF-3 and NFκB (NFκB p50 and NFκB p65) activity in the immune system. IKKε is a noncanonical kinase that phosphorylates inhibitors of NFκB, thus leading to dissociation and subsequent degradation of the inhibitor from the NFκB complex. IKKε is an oncogene that is often amplified in breast cancer. IKKε also functions to protect cells against apoptosis resulting from the DNA damage response.

**Protein Details:** Recombinant human IKKε protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids M1-V716, accession number NM\_014002 and N-terminally fused to GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

**Application Notes:** Recombinant IKKε is suitable for kinase assays and Western blot. The molecular weight of the protein is ~110.059 kDa. The activity of the protein is ~ 8 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: Casein 10 µg/50 µl, Recombinant IKKε : 100 ng/50 µl.

Kinase activity may vary depending on the substrate and reaction conditions used.

**Storage and Guarantee:** This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

