

## Data Sheet

### IKK $\beta$ , GST-tag

Human, recombinant, N-terminal GST-tag  
**Catalog#:** 40304  
**Lot#:** 150317-G1      **Conc.:** 0.37 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 3 mM DTT  
and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

#### References:

1. Woronicz, J.D. *et al.*, *Science*. 1997; **278**: 866-869.
2. Yin, M.J., *et al.*, *Nature*. 1998; **396**: 77-80.

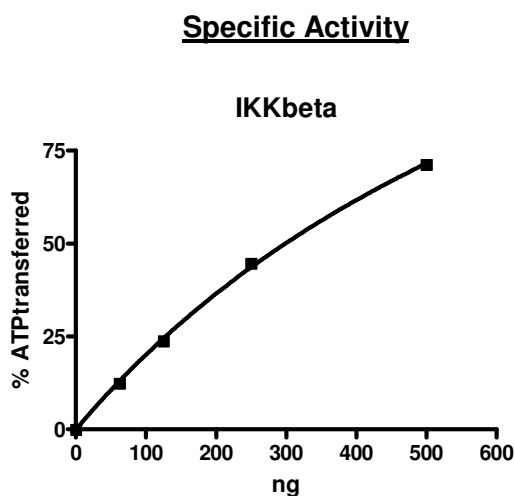
**Description:** Human IKK $\beta$ , also known as IKK2, IKKB, IKK $\beta$ , and NFKB1KB, GenBank Accession No. NM\_001556, full-length expressed in a baculovirus-infected Sf9 insect cell expression system, MW = 113 kDa.

**Specific Activity:**  $\geq 40$  pmol/min/ $\mu\text{g}$

**Assay Conditions:** Assay was done in a Kinase buffer using IKKtide as a substrate with 20  $\mu\text{M}$  ATP at  $30^{\circ}\text{C}$  for 45 min. The amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

**Application:** Useful for studying enzyme kinetics and screening small molecule inhibitors.

### Quality Assurance

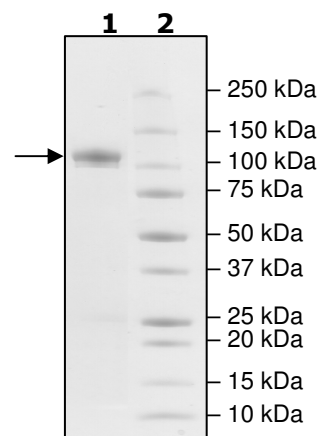


#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2  $\mu\text{g}$  IKKbeta

**Lane 2:**  
Protein Marker

**MW:** 113 kDa  
**Purity:**  $\geq 80\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)