

## Data Sheet

### ZAP70, Active

Human, recombinant, N-terminal GST tag

**Catalog #:** 40490

**Lot #:** 140813-1                      **Conc.:** 0.27 mg/ml

**Formulated in:** 40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% glycerol and 1.6 mM glutathione

**Stability:** >6 months at -80°C. *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

**References:**

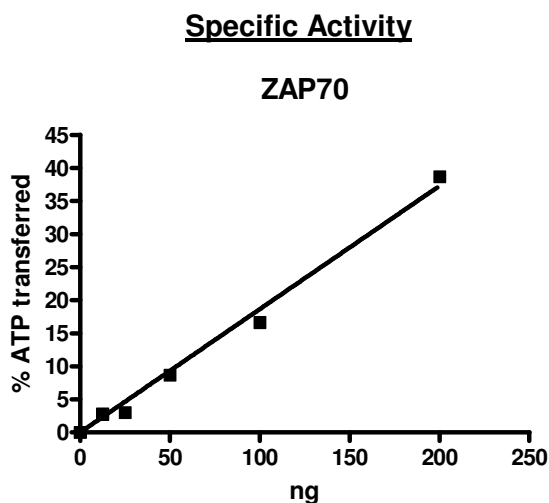
1. Ishaq, M., *et al. J Immunol.* 2007; **179(6)**:3851-8.
2. Siggs, O., *et al. Immunity.* 2007; **27(6)**:912-26.

**Description:** Human ZAP70, also known as ZAP-70, zeta-chain associated protein kinase, syk-related tyrosine kinase, SRK, or STD, GenBank Accession No. NM001079, full length with N-terminal GST tag, MW = 96 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 20 pmole/min/μg  
 Assay Conditions: Assay was done in Kinase buffer containing 0.2 mM DTT using Poly-(Glu4:Tyr) substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

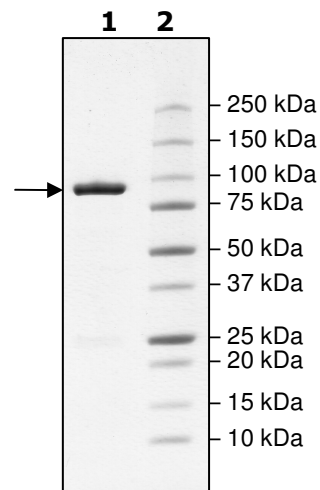
## Quality Assurance



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

**Lane 1:**  
 2.3 μg  
**Lane 2:**  
 Protein Marker

**MW:** 96 kDa  
**Purity:** ≥90%



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)