

## Data Sheet

### IRAK4, GST-Tag

Human, Recombinant, N-terminal GST tag  
**Catalog #:** 40064  
**Lot #:** 171113-3      **Conc.:** 1.21 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 300 mM NaCl, 2.4 mM KCl, 3 mM DTT, 10% Glycerol and 2 mM Glutathione.

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

**References:**

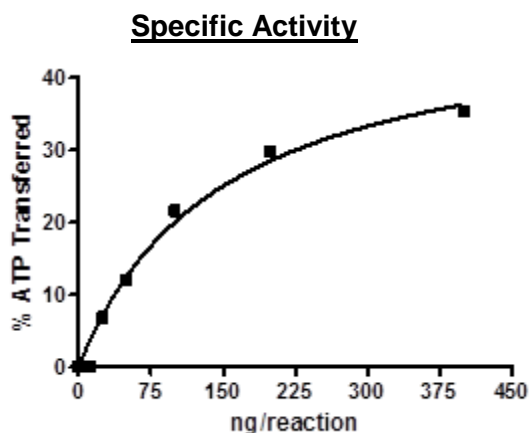
1. Ku C.L. *et al.*, *J. Exp. Med.* 2007; **204(10)**: 2407-2422.
2. Hoarau C. *et al.*, *J. Immunol.* 2007; **179(7)**: 4754-4765.

**Description:** Human interleukin-1 receptor-associated kinase 4 (IRAK4), GenBank Accession No. NM\_001114182, a.a 2-end, with N-terminal GST tag, MW = 78 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:**  $\geq 60$  pmol/min/ $\mu\text{g}$ .  
 Assay Conditions: Used 0.1 mg/ml MBP as a substrate and 20  $\mu\text{M}$  ATP. Reaction was done for 45 min at  $30^{\circ}\text{C}$ , and the amount of ATP was quantified using Kinase-Glo (Promega)

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

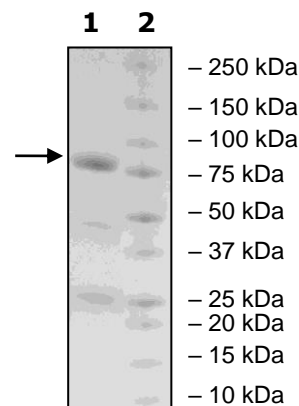
### Quality Assurance



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

**Lane 1:**  
2  $\mu\text{g}$  IRAK4  
**Lane 2:**  
Protein Marker

**MW:** 78 kDa  
**Purity:**  $\geq 67\%$



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 **1.858.259.5734**  
 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)