

## Data Sheet

### IRAK4, GST-Tag

Human, Recombinant, N-terminal GST-tag  
**Catalog #:** 40064  
**Lot #:** 171027-G      **Conc.:** 0.11 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 300 mM NaCl, 2.4 mM KCl, 3 mM DTT, 0.04% Tween-20 and 10% Glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**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.

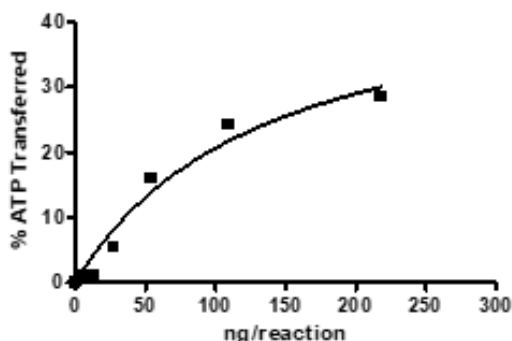
**Description:** Human interleukin-1 receptor-associated kinase 4 (IRAK4), GenBank Accession No. NM\_001114182, a.a 2-460(end), with N-terminal GST-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW = 78 kDa.

**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

### Specific Activity

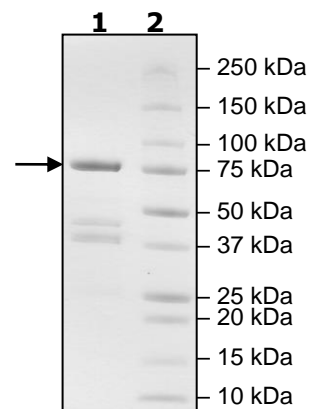


### 4-20% SDS-PAGE Coomassie staining

**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 **1.858.202.1401** 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)