

## Data Sheet

### IRAK1, His-tag

Human, Recombinant, N-terminal His-tag

**Catalog #:** 40202

**Lot #:** 180410-1

**Conc.:** 1.24 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 50 mM imidazole, and 20% glycerol.

**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.

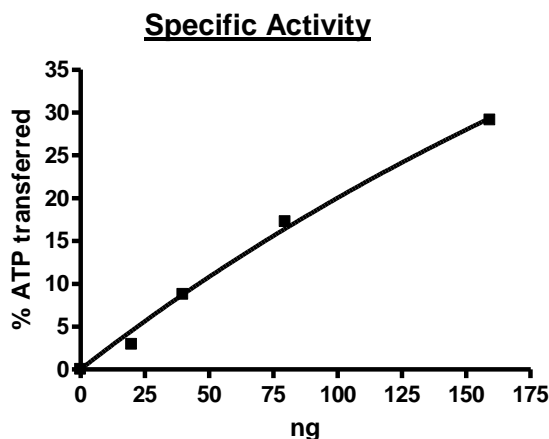
**Description:** Human interleukin-1 receptor-associated kinase 1 (IRAK1), GenBank Accession No. NM\_001569, a.a 186-530, with N-terminal His-tag, MW = 40 kDa, expressed in a Baculovirus infected Sf9 cell expression system.ss

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

**Assay Conditions:** Assay was done in Kinase buffer containing 1 mM DTT using MBP substrate (0.2 mg/ml) and 20  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  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.

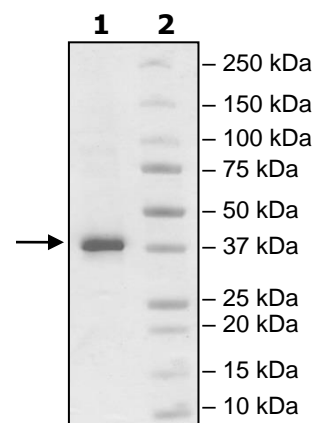
### Quality Assurance



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

**Lane 1:**  
2  $\mu\text{g}$  IRAK1  
**Lane 2:**  
Protein Marker

**MW:** 40 kDa  
**Purity:**  $\geq 90\%$



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 **1.858.259.5734**

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)