

Data Sheet

IRAK2

Human, recombinant, N-terminal GST tag

Catalog #: 40087

Lot#: 110520 **Conc.:** 0.1mg/ml

Formulated in: 50mM Tris-HCl, pH 7.5, 150 mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.1mM PMSF, 25% glycerol and 0.25 mM DTT.

Stability: >12months at -70°C

References:

1. Ku C.L. *et al.*, *J. Exp. Med.* **204 (10)**, 2407-2422 (2007).

2. Wan Y. *et al.*, *J Immunol.* 2011 Mar 1;**186(5)**:3006-14. Epub 2011 Jan 26.

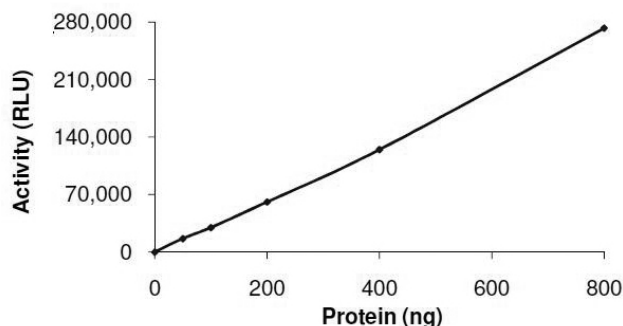
Description: Human IRAK2 (GenBank Accession No. NM_001570) full length with N-terminal GST tag, MW= 103 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 6 pmol/min/μg. Enzyme reaction is conducted in a buffer containing 50mM Tris-HCl, pH 7.5, 150 mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.1mM PMSF, 25% glycerol and 0.25 mM DTT.

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

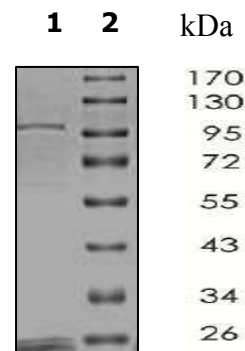
Quality Assurance

IRAK2 Specific Activity



10% SDS-PAGE
Coomassie staining

Lane 1: →
 IRAK2



MW: 103 kDa
Purity: ≥75%

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

Please visit our website at: www.bpsbioscience.com