

Product Information

IRAK1

 Product Number : **09-101**

Product description

Truncated human IRAK1 [194-712(end) amino acids of accession number NP_001560.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-IRAK1 was purified by using glutathione sepharose chromatography.

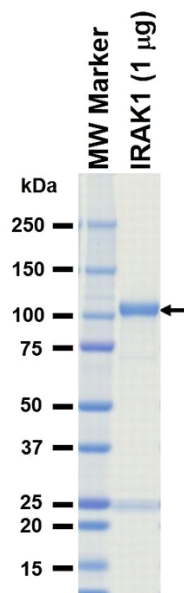
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
 Avoid repeating freeze-thaws.

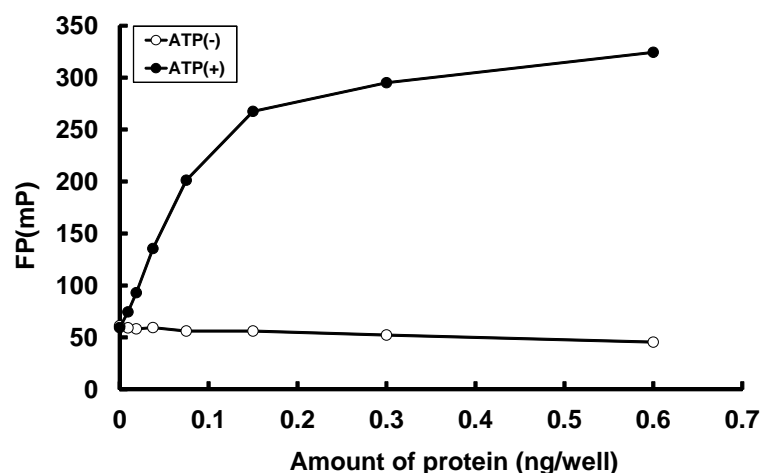
SDS-PAGE



Purity: 76 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : SRPKtide

ATP : 100 µM