

Data Sheet

BTK, GST-His-tag

Human, Recombinant, N-terminal GST-His tag

Catalog #: 40405

Lot #: 181019-GB **Conc.:** 0.05 mg/ml

Formulated in: 40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 KCl, 0.04% Tween-20 and 20% Glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

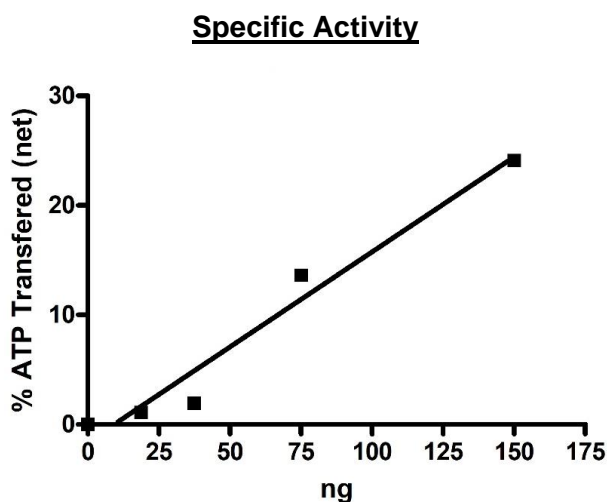
1. Liljeroos M., *et al.*, *Cell Signal*. 2007 Mar;**19(3)**:625-33
2. Jou ST., *et al.*, *Int. J. Immunogenet*. 2006 Oct;**33(5)**:361-369

Description: Human BTK also known as Bruton Tyrosine Kinase, ATK, BPK and Tyrosine-Protein Kinase BTK, GenBank Accession No. NM_000061, a.a 1-659(end) with N-terminal GST-His tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=106 kDa.

Specific Activity: 26 pmol/min/ μg .
 Assay Conditions: 40 mM Tris-HCl pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA and 1 mM DTT using 0.2 mg/ml Poly-(Glu4:Tyr)-biotin substrate and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent from Promega.

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

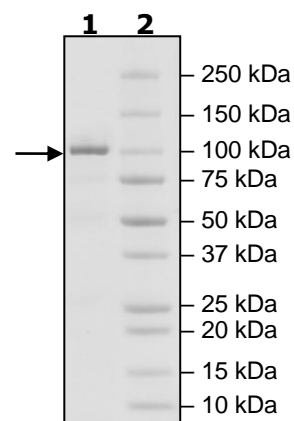
Quality Assurance



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

Lane 1:
2 μg BTK
Lane 2:
Protein Marker

MW: 106 kDa
Purity: $\geq 85\%$



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 **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com