



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

BTK

human, recombinant, N-terminal GST tag
Catalog #: 40405

Formulated in: 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

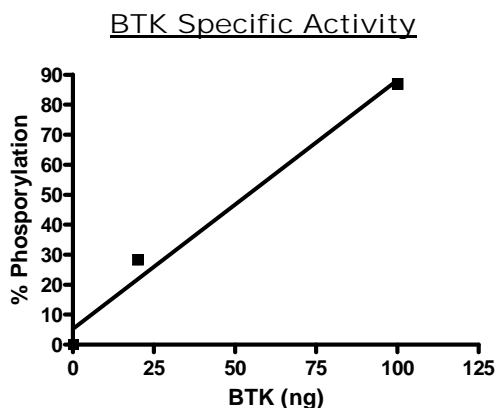
1. Liljeroos M., et al., Cell Signal. 2006 Aug 30
2. Jou ST., et al., Int J Immunogenet. 2006 Oct;33(5):361-369

Description: Human BTK (GenBank Accession No. NM_000061), full length with N-terminal GST tag, MW= 106 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

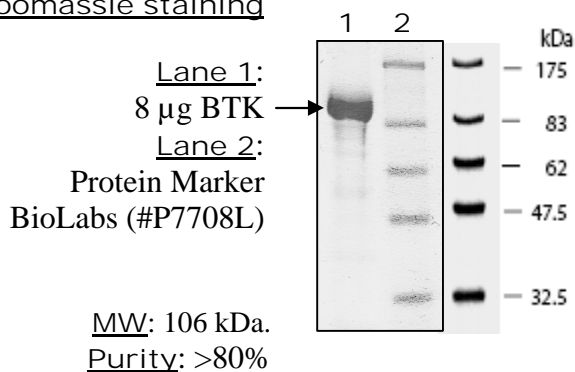
Specific Activity: 23 pmol/min/μg. Assay condition: The enzyme reaction was carried out for 1h at room temperature in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EDTA, 0.01% BRIJ-35 and 200 μM of ATP. Substrate : 2 μM Tyr peptide 1 from Invitrogen.

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

Quality Assurance



10% SDS-PAGE
Coomassie staining



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** or **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