

Data Sheet

DYRK1B, GST-tag

Human, Recombinant, N-terminal GST-tag
Catalog #: 100646
Lot #: 200317-G2 **Conc.:** 0.24 mg/ml

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

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

References:

1. Mao, J. *et al.*, *JBC*. **277(38)**: 35156-35161, 2002.
2. Guimera, J. W. *et al.*, *Genomics*. **57(3)**: 407-418, 1999.

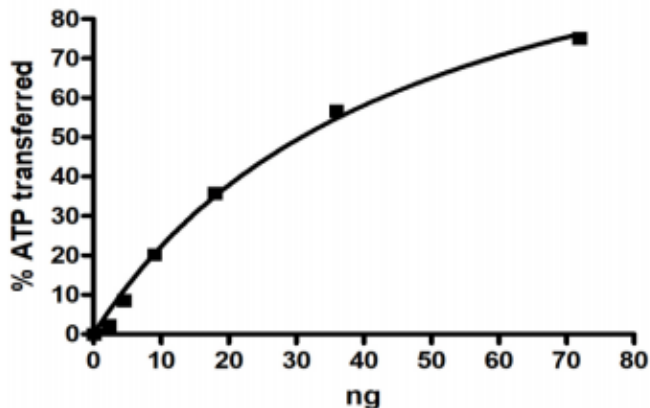
Description: Human Dual Specificity Tyrosine Phosphorylation Regulated Kinase 1B (DYRK1B), also known as Minibrain-Related Kinase (MIRK), GenBank Accession No. NM_004714, a.a. 2-629(end), with N-terminal GST-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW= 95 kDa.

Specific Activity: ≥500 pmole/min/μg
Assay Conditions: Assay was done in Kinase buffer containing DYRKtide substrate peptide 1 (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 35 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

Quality Assurance

Specific Activity

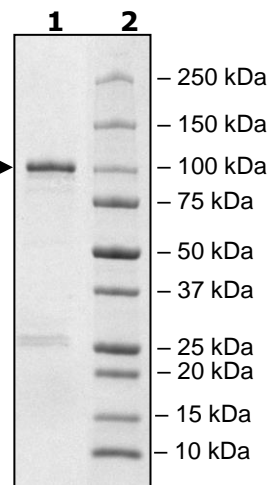


4-20% SDS-PAGE Coomassie staining

Lane 1:
2μg DYRK1B

Lane 2:
Protein Marker

MW: 95 kDa
Purity: ≥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