

Data Sheet

PYK2

Human, recombinant, N-terminal HIS tag

Catalog #: 40480

Lot #: 121203D **Conc.:** 0.26 mg/ml

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

Stability: >12 months at -80°C

References:

1. Cain R.J. et al., *J Cell Biol.* 2010 Mar 22;**188(6)**:863-76.
2. Collins M. et al., *Mol Immunol.* 2010 May;**47(9)**:1665-74.

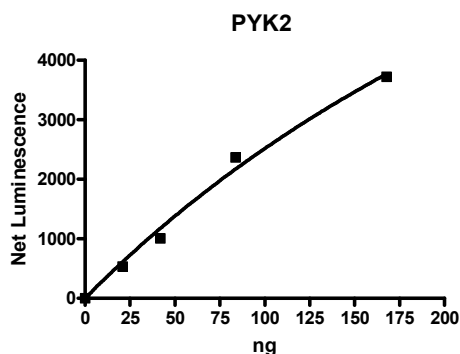
Description: Human PYK2 (GenBank Accession No. U33284), full length with N-terminal His tag, MW=120 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Specific Activity: 15 pmol/min/μg.
 Assay conditions: ADP-Glo assay with 0.2 mg/ml poly-(Glu4:Tyr) as a substrate and 100 μM ATP at final concentration. Reaction was done for 45 min at 30°C. ADP-Glo for 45 min at RT, and Kinase detection for 55 min at RT.

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

Quality Assurance

Specific Activity

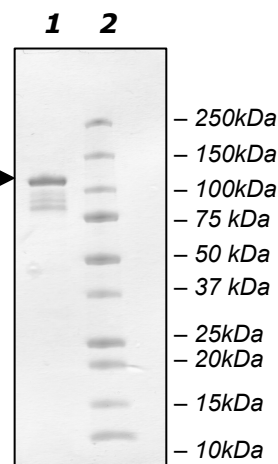


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
1.2 μg PYK2

Lane 2:
Protein Marker

MW: 120 kDa.
Purity: ≥62%



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 1.858.481.8694

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

Please visit our website at: www.bpsbioscience.com