

Data Sheet

PKLR Var 1 (PKR)

Human, recombinant, N-terminal His tag
 Catalog: 40500
 Lot#: 110201 Conc.: 0.76 mg/ml

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

Stability: >6 months at -80°C

References:

1. Bullido M.J. et al., *Neurobiol. Aging.*; **29(10):** 1160-6 (2008).
2. Russell S.T. et al., *Exp. Cell. Res.*; **315(1)** : 16-25 (2009).

Description:

PKR GenBank Accession No. NM_000298), full length with N-terminal His tag 2-574 (end), MW=63 kDa, expressed in an E. coli expression system.

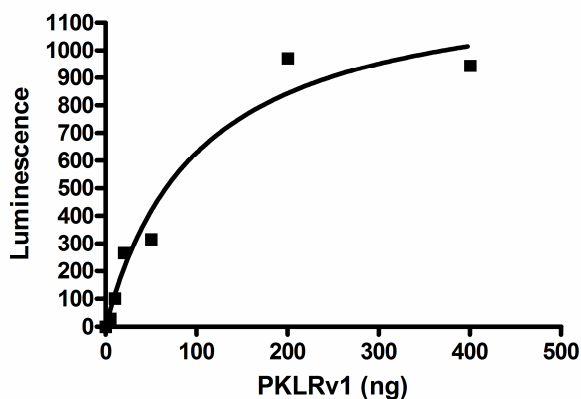
Specific Activity: 11 pmol/min/μg. Assay conditions: A 50 μl reaction is conducted in a buffer containing 50 mM Tris (pH 7.4), 10 mM MgCl₂, 50 mM KCl, 2 mM ADP, 10 mM phosphoenolpyruvate (PEP) at room temperature for 15 min. ATP production is detected using Kinase-Glo[®] Luminescent Kinase Assay Platform.

Application:

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

Quality Assurance

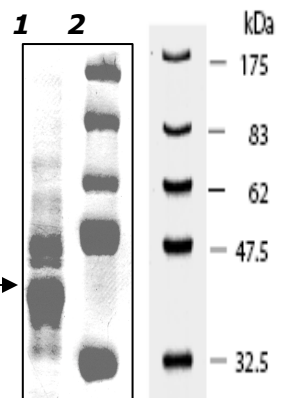
PKR Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
7.0 μg PKR
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 63 kDa
Purity: ≥43%



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