

Data Sheet

PKLR Var 1 (PKR)

Human, recombinant, N-terminal His tag
 Catalog: 40500
 Lot#: 110707G Conc.: 0.35 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.

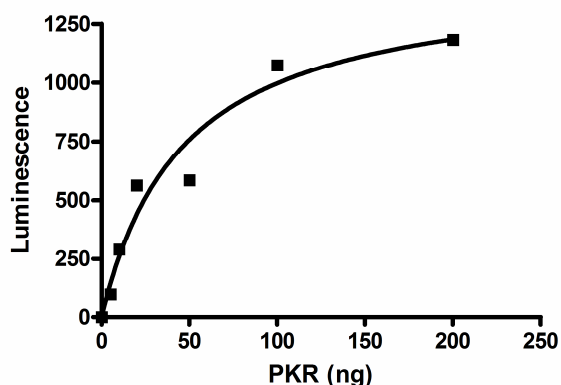
Specific Activity: 36 pmol/min/μg. Assay conditions: A 50 μl reaction is conducted in a buffer containing 50 mM Tris (pH 7.4), 10 mM MgCl₂, 50 mM KCl, 2 mM ADP, 10 mM phosphoenolpyruvate (PEP) at room temperature for 15 min. ATP production is detected using Kinase-Glo[®] Luminescent Kinase Assay Platform.

Application:

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

Quality Assurance

PKR Specific Activity

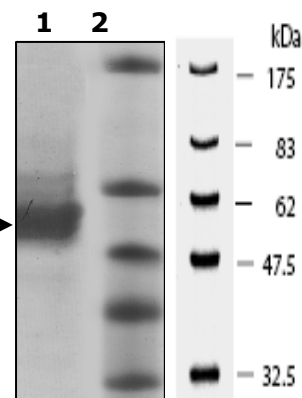


10% SDS-PAGE
Coomassie staining

Lane 1:
7.0 μg PKR

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 63 kDa
Purity: ≥80%



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