



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

BLK, Active

human, recombinant, N-terminal GST tag
Catalog #: 40401

Formulated in: 50mM Tris-HCl, pH 7.5,
150mM NaCl, 0.25mM DTT, 0.1mM EGTA,
0.1mM EDTA, 0.1mM PMSF, 25% glycerol.

Stability: >6 months at -80°C

References:

1. Dymecki, SM. et al. Science. 1990 Jan 19;247(4940):332-6.
2. Burkhardt, AL. et al. Proc Natl Acad Sci U S A. 1991 Aug 15;88(16):7410-4.

Description:

Recombinant full-length human BLK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is BC007371.

Specific Activity:

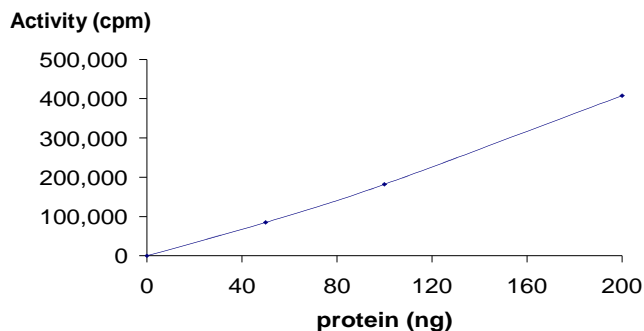
112 nmol/min/mg

Application:

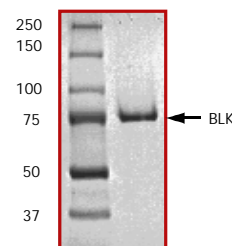
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

BLK Activity
112 nmol/min/mg



10% SDS-PAGE
Coomassie staining



MW: 84 kDa.
Purity: >90%

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