

Data Sheet

AKT2 Inactive

Human, recombinant, N-terminal His tag

Catalog #: 40001

Lot#: 100210 Conc.: 0.38 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 225 mM
imidazole, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Alessi *et al.*, *Current Biology*. **8: 69**, 1998
2. Cohen *et al.*, *FEBS Lett.* **410: 3**, 1997

Description: Human Akt2 (GenBank Accession No. NM_001626), full length with N-terminal HIS6 tag, MW=57 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

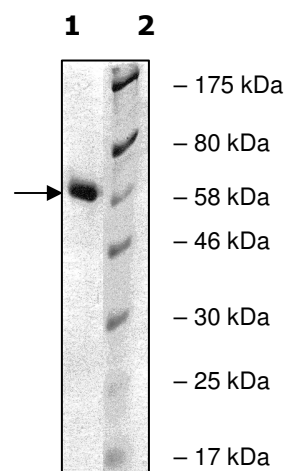
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
AKT2 Inactive
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 57 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com