



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

AKT2 Inactive

human, recombinant, , N-terminal His tag
Catalog #: 40001

Formulated in: 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

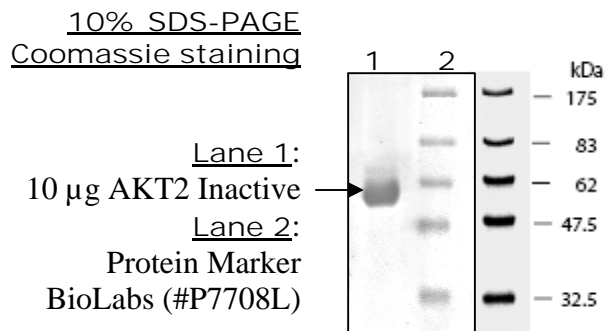
1. Alessi et al., Current Biology 8: 69, 1998
2. Cohen et al., FEBS Lett. 410: 3, 1997

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.

Specific Activity: Inactive

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

Quality Assurance



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 **1.858.202.1401** 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