



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

AKT1 Inactive

human, recombinant, , N-terminal His tag
Catalog #: 40000

Fomulated 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:

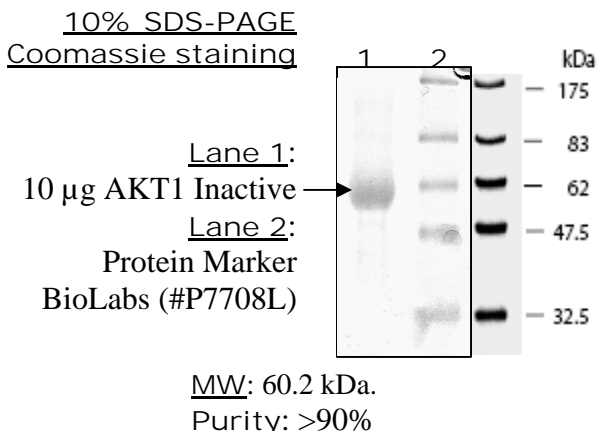
1. Alessi et al., Current Biology 8: 69, 1998
2. Cohen et al., FEBS Lett. 410: 3, 1997

Description: Human Akt1 (GenBank Accession No. NM_001014432), full length (a.a. 1-480) with N-terminal HIS6 tag, MW=60.2 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



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