

## Data Sheet

### AKT1 Inactive

Human, recombinant, N-terminal His tag

**Catalog #:** 40000

Lot#: 81222

Conc.: 1 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
124 mM NaCl, 2.2mM KCl, 10% glycerol,  
and 3 mM DTT.

**Stability:** >6 months at -80°C

**References:**

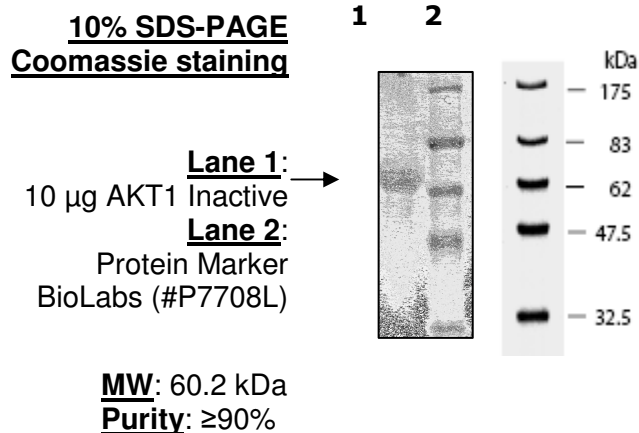
1. Alessi *et al.*, *Current Biology* **8: 69**, 1998
2. Cohen *et al.*, *FEBS Lett.* **410: 3**, 1997

**Description:** Human 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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)