



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### AKT2 active

Human, recombinant, N-terminal His tag

**Catalog #:** 40011

**Lot#** 110302-E1

**Conc.** 2 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 250 mM  
imidazole, 20% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

#### References:

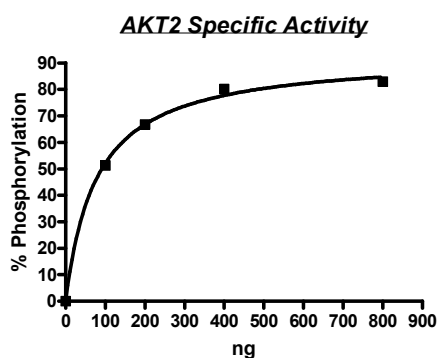
1. Alessi *et al.*, *Current Biology* **8**: 69, 1998
2. Cohen *et al.*, *FEBS Lett.* **410**: 3, 1997

**Description:** Recombinant full-length human AKT2 was co-expressed with PDK1 by baculovirus infection in Sf9 insect cells. The gene accession number is NM\_001626.

**Specific Activity:** 7 pmol/min/μg. The enzyme reaction is conducted in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl<sub>2</sub>, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35, and 2 μM substrate (SER/THR 6, Invitrogen) at 30 °C for 1 hour.

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

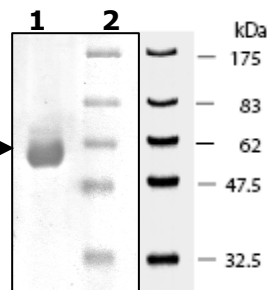
## Quality Assurance



### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
10 μg AKT2 active

**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](mailto:info@bpsbioscience.com)

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