

Data Sheet

AKT1, His-tag

Human, Recombinant, N-terminal His-tag

Catalog #: 40003

Lot#: 160504

Conc.: 0.27 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 200 mM Imidazole and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Kuijl C, *et al.*, *Nature*. 2007 Nov **29**;450 (7170):725-30.
2. Vogiatzi P and Giordano A. *Cancer Biol Ther*. 2007 Aug **3**;6(10).

Description:

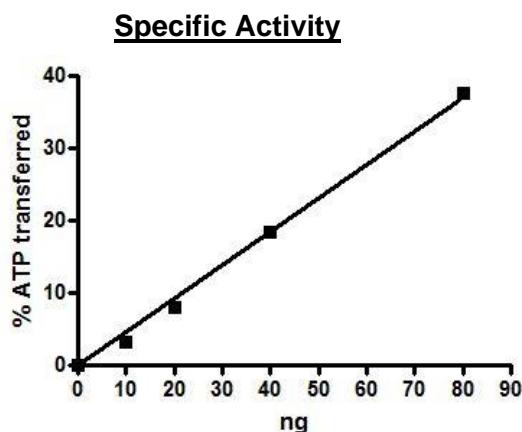
Human AKT Serine/Threonine Kinase 1 (AKT1), also known as Protein Kinase B Alpha (PKB α), and RAC, GenBank Accession No. NM_001014432, a.a. 1-480(full length) with N-terminal His-tag, was expressed in Sf9 insect cells via a baculovirus expression system. MW=60 kDa. The purified AKT1 is activated *in vitro* using PDK1 enzyme.

Specific Activity: >102 pmol/min/ μg .

Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using Crosstide substrate (0.1 mg/ml) and 20 mM ATP. Reaction was done at 30°C for 40 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

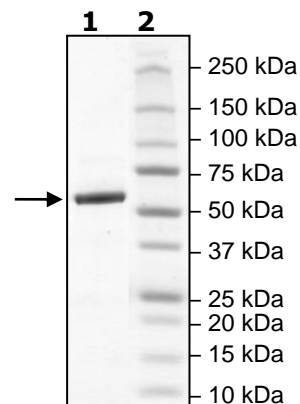
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg AKT1
Lane 2:
Protein Marker

MW: 60 kDa
Purity: $\geq 90\%$



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** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com