

Data Sheet

AKT3 active

human, recombinant, N-terminal His tag
Catalog #: 40012
Lot# 90212 **Conc.** 0.97 mg/ml

Fomulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2mM KCl, 200mM
 imidazole, 20% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

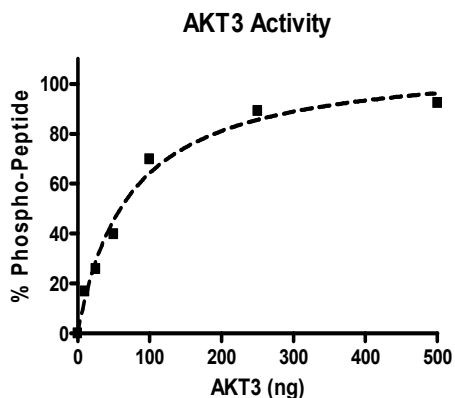
1. Coffey, P.J. et al. *Biochem J.* 1998 Oct 1; 335 (Pt 1):1-13.
2. Anderson, K.E. et al. *Curr Biol.* 1998 Jun 4;8(12): 684-91.

Description: Recombinant human AKT3(GenBank Accession No. NM_005465)(a-a2-end), with N-terminal HIS tag, MW= 60 kDa, expressed in baculovirus expression system.

Specific Activity: 14 pmol/min/ug. The enzyme reaction is conducted in a buffer containing 50 mM HEPES (pH7.5), 10mM MgCl₂, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35, and 2 μM substrate (SER/THR 6, Invitrogen) at room temperature for 1 hour.

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

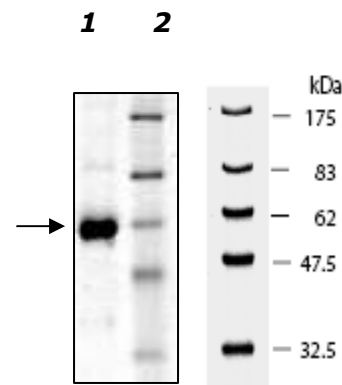
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 4.85 μg AKT3 active
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 60 kDa
Purity: ≥95%



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
 Please visit our website at: www.bpsbioscience.com