

# AKT2

aa107-481

**Product No.:** 0276-0000-2

**Lot:** 003

**Description:** Human AKT2

Amino acids A<sub>107</sub>-E<sub>481</sub> (as in GenBank entry NM\_001626.2)\*, N-terminally fused to GST-HIS<sub>6</sub>-Thrombin cleavage site

\*Sequence may contain documented polymorphisms  
Detailed sequence on request

**Product identity:** AKT2, Lot 003, was confirmed as AKT2 by specific Western Blotting using anti AKT2 antibody

**Theoretical MW**<sub>Fusion Protein</sub>: 72,844 Da

**Expression:** Baculovirus infected Sf9 cells

**Activation:** With PDK1  
(GenBank accession No.: NM\_002613)

**Purification:** One-step affinity purification using GSH-agarose

**Storage buffer:** 50 mM Tris-HCl, pH 8.0;  
100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.367 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Method for determination of K<sub>m</sub> value & Specific activity:**

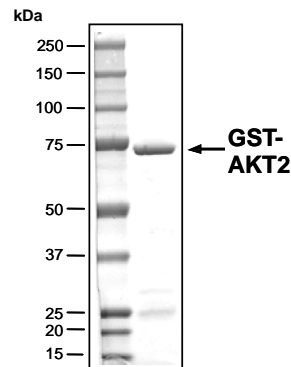
- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 3 µM Na-orthovanadate
  - 1.2 mM DTT
  - 2.5 µg / 50 µl PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: R<sub>11</sub>-GSK3(14-27)  
(R<sub>11</sub>-SGRARTSSFAEPGGK), 5 µg / 50 µl
  - Recombinant AKT2: 100 ng / 50 µl

- Filter binding assay  
MAPH membrane (Millipore)

**Specific activity:** 272 pmol/µg×min

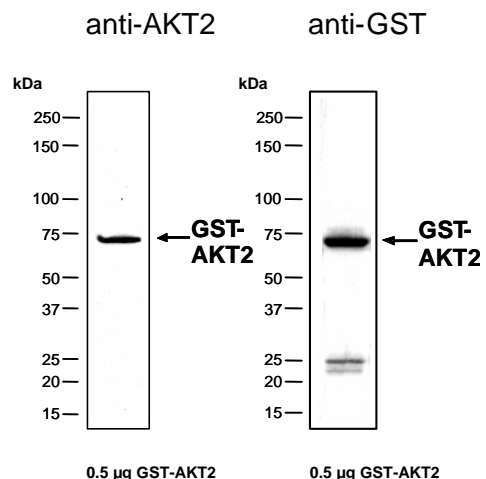
This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.

**Coomassie stain:**



2.0 µg GST-AKT2

**Western blot analysis:**



0.5 µg GST-AKT2

0.5 µg GST-AKT2

**Determination of K<sub>m</sub> value for ATP:**

