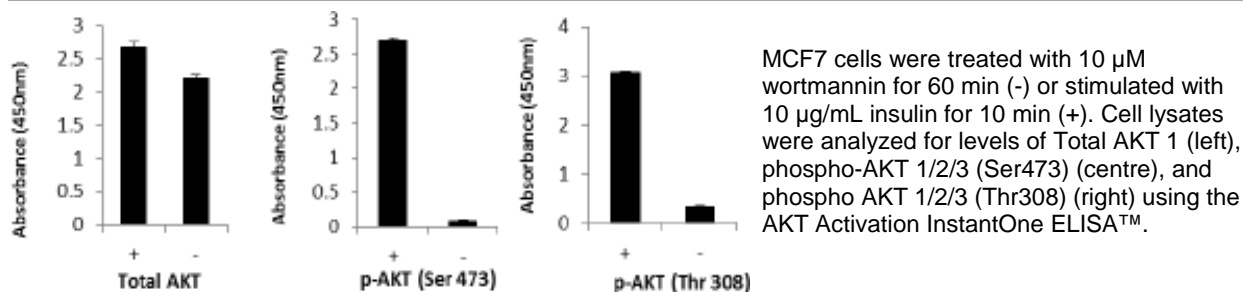


## AKT Activation InstantOne ELISA™ Kit

**Catalog Number:** 85-86049

**Also known as:** AKT, PKB, Protein kinase B, RAC-alpha serine/threonine-protein kinase

**RUO: For Research Use Only. Not for use in diagnostic procedures.**



### Product Information

**Contents:** AKT Activation InstantOne ELISA™

**Catalog Number:** 85-86049



**Temperature Limitation:** Store at 2-8°C.

**Batch Code:** Refer to vial

**Use By:** Refer to vial

### Description

The InstantOne ELISA™ is specifically engineered for accurate measurement of human, mouse, and hamster total AKT 1 and phosphorylated AKT 1/2/3 (Thr308 and Ser473) in cell lysates. Detection of phosphorylated rat AKT 1/2/3 (Thr308 and Ser473) and total rat AKT 1 is also expected. The InstantOne ELISA™ kit allows for fast analysis of samples in approximately one hour. All reagents used in a traditional sandwich ELISA are added in solution to a plate followed by a wash step and detection with the TMB colorimetric substrate.

AKT, which exists as multiple isoforms, is one of the principle kinases activated by phosphoinositide 3-kinase (PI3K). Activation of PI3K results in the generation of phosphatidylinositol (3,4,5)-triphosphate (PIP3). These lipid second messengers bind to the pleckstrin homology domain of AKT to promote its translocation to the plasma membrane for activation via phosphorylation at Thr308 and Ser473 by PDK1 and the mTOR TORC2 complex, respectively. Phosphorylation at both these sites is required for full activation of AKT Ser/Thr kinase activity. AKT phosphorylates over 50 known substrates, including GSK3, AS160, PRAS40, TSC 1, TCS 2, Raf-1, Bad, the FOXO family of transcription factors, and PFK2. Due to its established critical role in numerous biological processes, AKT is a frequent target for oncology-related drug development.

### Components

User Manual

AKT 1 (Total) Capture Antibody Reagent (Part No. IOAKTBCAR1T) - 1 mL

AKT 1 (Total) Detection Antibody Reagent (Part No. IOAKTBDAR1T) - 1 mL

phospho-AKT 1/2/3 (Thr308) Capture Antibody Reagent (Part No. IOPAKTBCAR2T) - 1 mL

phospho-AKT 1/2/3 (Thr308) Detection Antibody Reagent (Part No. IOPAKTBDAR2T) - 1 mL

phospho-AKT 1/2/3 (Ser473) Capture Antibody Reagent (Part No. IOPAKTACAR1T) - 1 mL

phospho-AKT 1/2/3 (Ser473) Detection Antibody Reagent (Part No. IOPAKTADAR1T) - 1 mL

InstantOne™ Assay Plate (Part No. IOAP96) - 96 well (sealed)

Cell Lysis Buffer (5X) (Part No. IOCLB1) - 10 mL

Enhancer Solution (Part No. IOES1) - 1 mL

Wash Buffer (10X) (Part No. IOWB1) - 15 mL

Detection Reagent (TMB Substrate) (Part No. IODR1) - 12 mL

Stop Solution (Part No. IOSS1) - 12 mL

Positive Control Cell Lysate (lyophilized) (Part No. IOAKTAPC1) - 0.25 mL

Plate seal - 1 ea

Not for further distribution without written consent. Copyright © 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Tel: 1 800 9556288 • [thermofisher.com/askaquestion](http://thermofisher.com/askaquestion) • [thermofisher.com/support](http://thermofisher.com/support)

---

## AKT Activation InstantOne ELISA™ Kit

**Catalog Number:** 85-86049

**Also known as:** AKT, PKB, Protein kinase B, RAC-alpha serine/threonine-protein kinase

**RUO: For Research Use Only. Not for use in diagnostic procedures.**

---

### Applications Reported

This ELISA antibody has been reported for use in ELISA.

### Applications Tested

This assay has been validated for the detection of total AKT 1 and phosphorylated AKT 1/2/3 (Thr308 and Ser473) in stimulated MCF7 cell lysates (data shown above).

### Special Notes

Reagents for 32 wells of total AKT1, 32 wells of phospho-AKT1/2/3 (Thr308), and 32 wells of phospho-AKT1/2/3 (Ser473).

### Related Products

85-86042 phospho-AKT 1/2/3 (Ser473) InstantOne ELISA™

85-86044 phospho-AKT 1/2/3 (Thr308) InstantOne ELISA™

85-86046 AKT (Total/Phospho) InstantOne ELISA™

85-86047 AKT 1 (Total) InstantOne ELISA™

IOCLB InstantOne ELISA Cell Lysis Buffer (5X) – 100 mL

IOES InstantOne ELISA Enhancer Solution – 100 mL

IOWB InstantOne ELISA Wash Buffer (10X) – 500 mL