

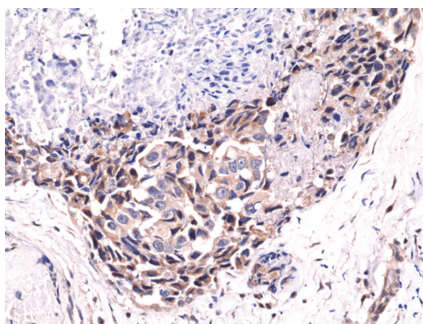
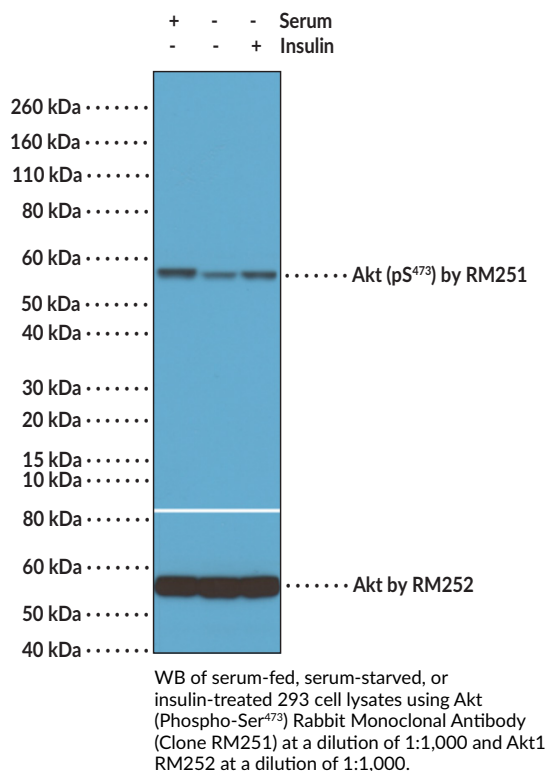
PRODUCT INFORMATION

Akt (Phospho-Ser⁴⁷³) Rabbit Monoclonal Antibody (Clone RM251) Item No. 42814

Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	PKB, Protein Kinase B, RAC Serine/threonine-protein Kinase, RAC-PK
Immunogen:	A phosphopeptide corresponding to human Akt (phospho-Ser ⁴⁷³)
Cross Reactivity:	(+) Akt1/2/3 (phospho-Ser ⁴⁷³); (-) Unmodified Akt
Species Reactivity:	(+) Human, mouse, rat, bovine
Molecular Weight:	56 kDa
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM251
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:200-1:500 and 1-1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue sections using Akt (Phospho-Ser⁴⁷³) Rabbit Monoclonal Antibody (Clone RM251) at a dilution of 1:400.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/12/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Akt, also known as protein kinase B (PKB), is a serine/threonine kinase belonging to the AGC kinase family and has three isoforms in mammals: Akt1, Akt2, and Akt3.^{1,2} Akt kinases function downstream of activated tyrosine kinases and PI3K to regulate a variety of cellular processes, including cell size, growth, proliferation, and survival, as well as genome stability, glucose metabolism, and neovascularization.² Akt is composed of an N-terminal pleckstrin homology (PH) domain, which binds to phosphatidylinositol-(3,4,5)-triphosphate (PIP₃) and phosphatidylinositol-(3,4)-diphosphate (PIP₂), a kinase domain, and a C-terminal regulatory hydrophobic motif. Akt1 and Akt2 are ubiquitously expressed whereas Akt3 is primarily expressed in the placenta, ovary, and spleen, and each isoform shows distinct subcellular localizations: Akt1 in the cytoplasm, Akt2 in the mitochondria, and Akt3 in the nucleus and nuclear membrane.^{3,4} Akt is activated following phosphorylation at threonine 308, 309, and 305 and serine 473, 474, and 472 for Akt1, -2, and -3, respectively.⁴ Hyperactivation of Akt *via* mutations, amplification, or overexpression is associated with various cancers.² Cayman's Akt (Phospho-Ser⁴⁷³) Rabbit Monoclonal Antibody (Clone RM251) can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes Akt (phospho-Ser⁴⁷³) at 56 kDa from human, mouse, rat, and bovine samples.

References

1. Dummer, B. and Hemmings, B.A. Physiological roles of PKB/Akt isoforms in development and disease. *Biochem. Soc. Trans.* **35**(Pt 2), 231-235 (2007).
2. Bellacosa, A., Kumar, C.C., Di Cristofano, A., *et al.* Activation of AKT kinases in cancer: Implications for therapeutic targeting. *Adv. Cancer Res.* **94**, 29-86 (2005).
3. Masure, S., Haefner, B., Wesselink, J.J., *et al.* Molecular cloning, expression and characterization of the human serine/threonine kinase Akt-3. *Eur. J. Biochem.* **265**(1), 353-360 (1999).
4. Santi, S.A. and Lee, H. The Akt isoforms are present at distinct subcellular locations. *Am. J. Physiol. Cell Physiol.* **298**(3), C580-C591 (2010).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM