

Product Information

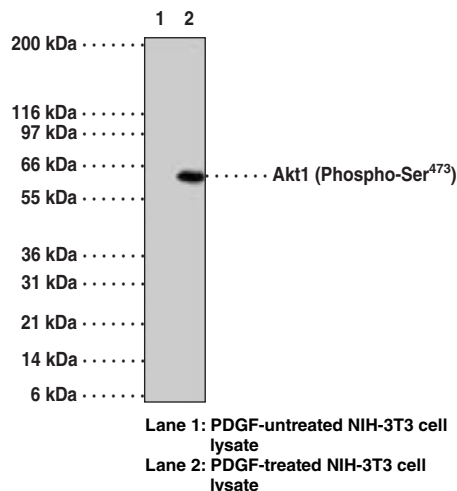


Akt1 (Phospho-Ser473) Monoclonal Antibody (Clone 104A282)

Item No. 13733

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Antigen:** Synthetic peptide from human Akt1 containing phospho-serine⁴⁷³
- Host:** Mouse, clone 104A282
- Isotype:** IgG_{1κ}
- Cross Reactivity:** (+) Human and mouse Akt1; predicted to cross react with chicken and frog Akt1
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Application:** Immunoprecipitation (IP) and western blot (WB). The recommended starting concentration for WB is 2 µg/ml and 1 µg/ml for IP.

Akt, protein kinase B (PKB), is a serine/threonine kinase which is involved in many cellular signaling pathways and acts as a transducer of many functions initiated by growth factor receptors that activate phosphatidylinositol 3-kinase (PI3-kinase). The major activity of Akt/PKB is to mediate cell survival. Akt/PKB is also believed to be a critical factor in the genesis of cancer, as the tumor suppressor PTEN was found to antagonise PI3 kinase and Akt/PKB kinase activity. Akt/PKB phosphorylation is critical for its activity. The major phosphorylation sites required for Akt activation have been identified as threonine 308 and serine 473. Serine⁴⁷³ is phosphorylated by MAPKAP kinase 2.



Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13733

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 04/10/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com