

Product Information



Ribosomal S6 Kinase 2 Polyclonal Antibody

Catalog No. 10009411

Synonym:	RSK2
Supplied as:	100 µl of affinity-purified antibody in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg BSA per ml and 50% glycerol.
Host:	Rabbit
Antigen:	Peptide corresponding to amino acid residues from the C-terminal region of rat ribosomal S6 kinase 2.
Cross Reactivity:	(+) Rat RSK2; expected to react with bovine, canine, chicken, murine, and human RSK2
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

The p90 ribosomal S6 kinases (RSK) 1-4 are downstream members of the extracellular signal-regulated kinase (ERK)/MAPK cascade. The loss of RSK2 activity in humans leads to Coffin-Lowry syndrome, which is characterized by mental retardation and growth deficits.¹ Recent work suggests that RSK2 exerts a tonic regulation on G protein-coupled signalling.²

References

1. Hanauer, A. and Young, I.D. Coffin-Lowry syndrome: Clinical and molecular features. *J. Med. Genet.* **39**, 705-713 (2002).
2. Sheffler, D.J., Kroeze, W.K., Garcia, B.G., *et al.* p90 ribosomal S6 kinase 2 exerts a tonic brake on G protein-coupled receptor signaling. *Proc. Natl. Acad. Sci. USA* **103**, 4717-4722 (2006).

Related Products

p38 MAPK (Phospho-Thr¹⁸⁰/Tyr¹⁸²) Polyclonal Antibody - Cat. No. 10009177 • MEK 1/2 (Phospho-Ser^{218,222}) Polyclonal Antibody - Cat. No. 10009178 • ERK/MAPK (Phospho-Thr²⁰²/Tyr²⁰⁴) Polyclonal Antibody - Cat. No. 10009179 • GSK3β (Phospho-Ser⁹) Polyclonal Antibody - Cat. No. 10009374

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

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

MATERIAL 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 Material 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, 01/05/2010