

PRODUCT INFORMATION



Ribosomal S6 Kinase 2 Polyclonal Antibody

Item No. 10009411

Overview and Properties

Contents:	This vial contains 100 µl of peptide affinity-purified antibody.
Synonym:	RSK2
Immunogen:	Peptide from the C-terminal region of rat RSK2
Species Reactivity:	(+) Rat RSK2; expected to react with bovine, canine, chicken, mouse, and human RSK2
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly
Storage Buffer:	10 mM HEPES, pH 7.5, containing 150 nM sodium chloride, 100 µg BSA/ml and 50% glycerol
Host:	Rabbit
Application:	Western blot; the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

The p90 ribosomal S6 kinases (RSKs) 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 signaling.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).

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, 08/11/2016

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