

PRODUCT INFORMATION



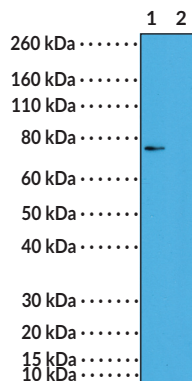
RSK1 (Phospho-Thr³⁵⁹/Ser³⁶³) Rabbit Monoclonal Antibody (Clone RM233)

Item No. 32193

Overview and Properties

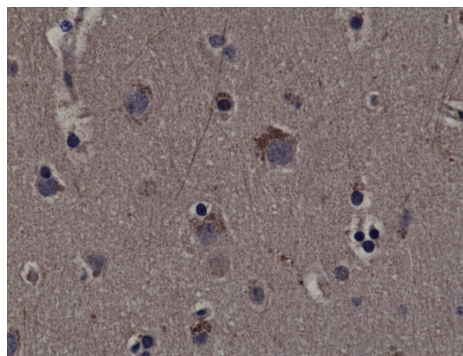
Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonym:	p90 Ribosomal S6 Kinase 1 (phospho-T ³⁵⁹ /S ³⁶³)
Immunogen:	Peptide corresponding to RSK1 (phospho-Thr ³⁵⁹ /Ser ³⁶³)
Cross Reactivity:	(+) RSK1 (phospho-Thr ³⁵⁹ /Ser ³⁶³); (-) RSK1 without phosphorylation at Thr ³⁵⁹ /Ser ³⁶³
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM233
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200-1:500 and 1:1,000-1:2,000 for IHC and WB, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: A431 cells treated
Lane 2: A431 cells untreated

WB of A431 cells treated with EGF or left untreated using RSK1 (Phospho-Thr³⁵⁹/Ser³⁶³) Rabbit Monoclonal Antibody (Clone RM233) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human cerebral cortex tissue using RSK1 (Phospho-Thr³⁵⁹/Ser³⁶³) Rabbit Monoclonal Antibody (Clone RM233) at a dilution of 1:200.

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, 01/19/2024

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

p90 ribosomal S6 kinase 1 (RSK1) is a member of the RSK family of serine/threonine kinases that mediates RAS/RAF/MEK/ERK/MAPK signaling.¹ It contains N- and C-terminal kinase domains connected by a linker region, which contains various residues, including threonine 359 (Thr³⁵⁹) and serine 363 (Ser³⁶³), that are subject to phosphorylation and lead to its activation. RSK1 is ubiquitously expressed and localized to the cytoplasm, where it is in a complex with its activator ERK1/2 in quiescent cells.^{1,2} Upon stimulation of receptor tyrosine kinases (RTKs) by a variety of cytokines, neurotransmitters, or hormones, RSK1 accumulates at the plasma membrane and is phosphorylated by ERK1/2 and 3-phosphoinositide-dependent protein kinase 1 (PDK1), resulting in dissociation of the ERK1/2-RSK1 complex and activation of RSK1.¹⁻³ RSK1 phosphorylates a variety of substrates that have roles in several cellular processes, including cell growth, survival, and proliferation.² Tumor RSK1 levels are increased in patients with breast or prostate cancer.¹ Cayman's RSK1 (Phospho-Thr³⁵⁹/Ser³⁶³) Rabbit Monoclonal Antibody (Clone RM233) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Casalvieri, K.A., Matheson, C.J., Backos, D.S., *et al.* Selective targeting of RSK isoforms in cancer. *Trends Cancer* **3(4)**, 302-312 (2017).
2. Romeo, Y., Zhang, X., and Roux, P.P. Regulation and function of the RSK family of protein kinases. *Biochem. J.* **441(2)**, 553-569 (2012).
3. Anjum, R. and Blenis, J. The RSK family of kinases: Emerging roles in cellular signalling. *Nat. Rev. Mol. Cell Biol.* **9(10)**, 747-758 (2008).

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