

Product Information



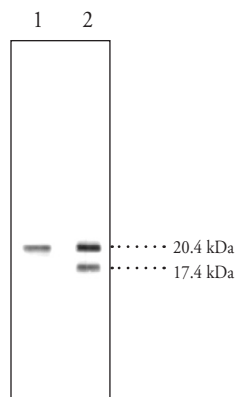
Caveolin 1/3 Polyclonal Antiserum

Catalog No. 100830

Contents:	100 µl of lyophilized antiserum
Host:	Rabbit
Stability:	≥1 year at -20°C
Antigen:	Rat caveolin-3 amino acids 19-41 (CKEIDLVNRPKNINEDIVKVDF) conjugated to KLH. This sequence is identical to human and murine caveolin-3 and shares significant identity with caveolin-1 of several species. Murine, rat, or human caveolin-3 aa 19-41 CKEIDLVNRPKNINEDIVKVDF Murine, rat, or human caveolin-1 aa 46-68 t KEIDLVNRPK h I N d D v V K i D F Zebrafish caveolin-1a aa 50-71 - KEIDLVNRPK h I N d D v V K V D F Zebrafish caveolin-1b aa 17-3 - KEIDLVNRPK h I N d D v V K V D F
Cross-reactivity:	(+) Caveolin-1 and caveolin-3 from most mammalian species. Also, it has not been tested but should detect caveolin-1 isoforms in some frogs (<i>X. laevis</i>) and fish like zebrafish (<i>D. rerio</i>) and pufferfish (<i>T. rubripes</i>)
Application:	Western blot; other applications not tested

Caveolin is an integral membrane protein and is a principle component of caveolea, flask-shaped plasma membrane invaginations.

Currently, 3 isoforms of caveolin (caveolin-1, caveolin-2, and caveolin-3), which differ in their primary sequence and tissue distribution, have been cloned.^{1,2,3} Caveolin-1, cloned from a variety of species, has 178 amino acids with a molecular mass of approximately 21 kDa. Caveolin-3 (rat) contains 151 amino acids with a predicted molecular mass of 17,404 Da.³ Based on protein sequence homology, rat caveolin-1 and caveolin-3 are approximately 65% identical. Caveolin-1 is found in adipocytes, endothelial cells, and fibroblasts, whereas Caveolin-3 is found predominately in muscle tissue.³



1. Bovine aorta endothelial cell lysate
2. Adult rat ventricular myocyte cell lysate

References

1. Tang, Z.L., Scherer, P.E., Lisanti, M.P. The primary sequence of murine caveolin reveals a conserved consensus site for phosphorylation by protein kinase C. *Gene* **147**, 299-300 (1994).
2. Scherer, P.E., Okamoto, T., Nishimoto, I., *et al.* Identification, sequence, and expression of caveolin-2 defines a caveolin gene family. *Proc. Natl. Acad. Sci. USA* **93**, 131-135 (1996).
3. Tang, Z.L., Scherer, P.E., Okamoto, T., *et al.* Molecular cloning of caveolin-3, a novel member of the caveolin gene family expressed predominantly in muscle. *J. Biol. Chem.* **271**, 2255-2261 (1996).

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSDS supervisor at 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, 05/11/2009

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