

# PRODUCT INFORMATION



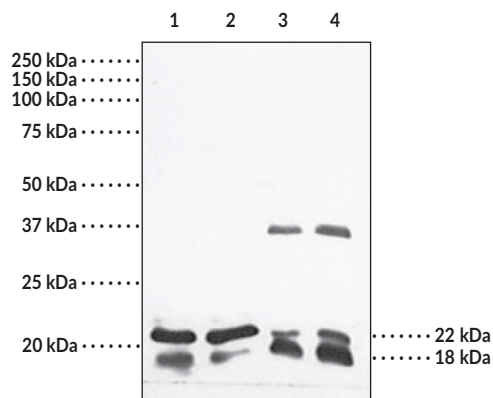
## Caveolin-1/3 Polyclonal Antiserum

Item No. 100830

### Overview and Properties

<b>Contents:</b>	This vial contains polyclonal antiserum
<b>Synonym:</b>	Cav-1/3
<b>Immunogen:</b>	Synthetic peptide from the N-terminal region of rat caveolin-3
<b>Cross Reactivity:</b>	(+) Caveolin-1 and caveolin-3
<b>Species Reactivity:</b>	(+) Human, murine, and rat; other species not tested
<b>Uniprot No.:</b>	P51638
<b>Form:</b>	Lyophilized
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	Polyclonal antiserum when reconstituted in 100 µl of double distilled water
<b>Host:</b>	Rabbit
<b>Applications:</b>	Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: Mouse skeletal muscle supernatant (30 µg)  
Lane 2: Mouse skeletal muscle supernatant (50 µg)  
Lane 3: Mouse heart homogenate (30 µg)  
Lane 4: Mouse heart homogenate (50 µg)

#### 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, 11/02/2023

#### 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

---

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.<sup>1,2,3</sup> 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.<sup>3</sup> 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.<sup>3</sup>

## 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(2)**, 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(1)**, 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(4)**, 2255-2261 (1996).

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