

PRODUCT INFORMATION



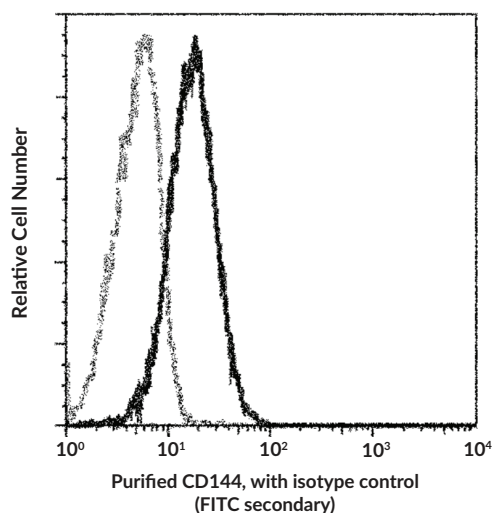
VE-Cadherin/CD144 Monoclonal Antibody (Clone 01)

Item No. 37086

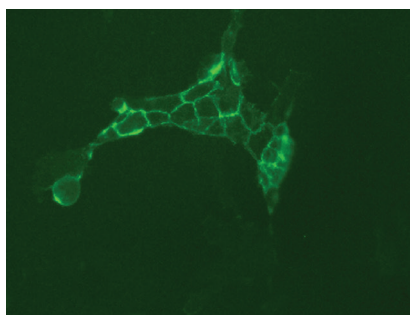
Overview and Properties

Contents:	This vial contains 50, 100, or 200 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	7B4 Antigen, Cadherin-5, Vascular Endothelial Cadherin
Immunogen:	Recombinant human VE-cadherin
Cross Reactivity:	(+) VE-cadherin
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	01
Host:	Mouse
Isotype:	IgG1
Applications:	Flow cytometry (FC), Immunocytochemistry (ICC), and Immunofluorescence (IF); the recommended starting dilution for FC is 1:25-1:100 and 1:10-1:50 for ICC and IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Flow cytometric analysis of human VE-cadherin/CD144 expression on HUVEC cells. Cells were stained with purified VE-Cadherin/CD144 Monoclonal Antibody (Clone 01) then a FITC-conjugated secondary antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of VE-cadherin in HUVEC cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with VE-Cadherin/CD144 Monoclonal Antibody (Clone 01) (dilution ratio 1:30) at 4°C overnight. Then cells were stained with an Alexa Fluor®488-conjugated goat anti-mouse IgG secondary antibody (green). Positive staining was localized to the cell membrane.

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, 02/13/2024

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

VE-cadherin, also known as CD144, is a type II classical cadherin and member of the cadherin superfamily of calcium-dependent cell adhesion molecules.¹ It is composed of an N-terminal signal peptide, a pro-peptide, five extracellular cadherin repeats that bind to homotypic cadherins, a transmembrane domain, and a cytoplasmic tail that interacts with the cytoskeleton. VE-cadherin is endothelial cell-specific and is localized to intracellular junctions where it participates in vascular structure assembly.²⁻⁴ An anti-VE-cadherin antibody in combination with an anti-CD31 antibody inhibits tube formation in human umbilical vein endothelial cells (HUVECs).⁵ Knockout of *Cdh5*, the gene encoding VE-cadherin, in mouse embryonic stem cells prevents the formation of vessel-like structures in 11-day-old embryoid bodies.³ VE-cadherin is overexpressed in melanoma and breast cancer and is positively correlated with poor prognosis.⁶ Cayman's VE-Cadherin/CD144 Monoclonal Antibody (Clone 01) can be used for flow cytometry (FC), immunocytochemistry (ICC), and immunofluorescence (IF) applications.

References

1. Yu, W., Yang, L., Li, T., *et al.* Cadherin signaling in cancer: Its functions and role as a therapeutic target. *Front Oncol.* **9**, 989 (2019).
2. Lampugnani, M.G., Resnati, M., Raiteri, M., *et al.* A novel endothelial-specific membrane protein is a marker of cell-cell contacts. *J. Cell Biol.* **118(6)**, 1511-1522 (1992).
3. Vittet, D., Buchou, T., Schweitzer, A., *et al.* Targeted null-mutation in the *vascular endothelial-cadherin* gene impairs the organization of vascular-like structures in embryoid bodies. *Proc. Natl. Acad. Sci. USA* **94(12)**, 6273-6278 (1997).
4. Breviario, F., Caveda, L., Corada, M., *et al.* Functional properties of human vascular endothelial cadherin (7B4/cadherin-5), an endothelium-specific cadherin. *Arterioscler. Thromb. Vasc. Biol.* **15(8)**, 1229-1239 (1995).
5. Matsumura, T., Wolff, K., and Petzelbauer, P. Endothelial cell tube formation depends on cadherin 5 and CD31 interactions with filamentous actin. *J. Immunol.* **158(7)**, 3408-3416 (1997).
6. Bartolomé, R.A., Torres, S., Isern de Val, S., *et al.* VE-cadherin RGD motifs promote metastasis and constitute a potential therapeutic target in melanoma and breast cancers. *Oncotarget* **8(1)**, 215-227 (2017).

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