

PRODUCT INFORMATION

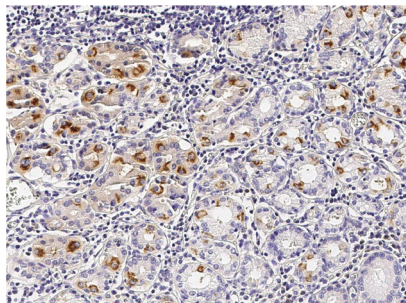
VE-Cadherin/CD144 Rabbit Monoclonal Antibody (Clone 012)

Item No. 37085

Overview and Properties

Contents:	This vial contains 50 or 100 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; other species not tested
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	012
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC); the recommended starting concentration is 1:250-1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemical staining of formalin-fixed and paraffin-embedded human stomach using VE-Cadherin/CD144 Rabbit Monoclonal Antibody (Clone 012) at a dilution of 1:500.

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 Rabbit Monoclonal Antibody (Clone 012) can be used for immunohistochemistry (IHC).

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**(16), 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