

# VE-cadherin (human) monoclonal antibody (BV6)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 18

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-803-3050	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Vascular endothelial cadherin, Cadherin-5, CD144, 7B4 antigen
<b>Application</b>	ELISA, Flow Cytometry, ICC, IHC (FS), IP, WB
<b>Application Notes</b>	Detects a band of ~130-140kDa by Western blot.
<b>Clone</b>	BV6
<b>Crossreactivity</b>	Does not cross-react with other species.
<b>Formulation</b>	Liquid. In PBS containing 1mg/ml BSA and 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Human endothelial cells.
<b>Isotype</b>	IgG2a
<b>Purity Detail</b>	Affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Immunocytochemistry (fluorescence and peroxidase staining)Western Blot (non-reducing)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human

UniProt ID

P33151

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)