

# CD54 (human) (domain D1) monoclonal antibody (RR1/1)

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

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-805-016-C050	Purified - 50µg
------------------	-----------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ICAM-1, Intercellular adhesion molecule-1
Application	ELISA, Flow Cytometry, FUNC, IHC (FS), WB
Application Notes	Detects a band of ~90kDa by Western blot. Functional Application: blocks adhesion of leukocytes to endothelial cells.
Clone	RR1/1
Formulation	Liquid. In PBS, pH 7.2-7.4.
Host	Mouse
Immunogen	Human CD54 (ICAM-1).
Isotype	IgG1
Recommendation	Flow Cytometry (suitable as primary antibody in staining for FACS
Dilutions/Conditions	analysis)Immunohistochemistry (stains acetone-fixed cryostat sections or cell smears. Alkaline phosphatase-anti-alkaline phosphatase (APAAP) procedure, peroxidase anti-peroxidase (PAP) procedure or the three stage immunoperoxidase technique are recommended; use at a concentration of 0.5-1.25µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes the D1 domain of CD54.



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

P05362

European Sales Office  
ENZO LIFE SCIENCES  
(EU) 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)

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee**.

For local distributors and detailed product information visit us online: [www.enzo.com](http://www.enzo.com) All product names, logos, and brands are the property of their respective owners.