

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

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

Citations: 2

[View Online »](#)

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 Dilutions/Conditions	Flow Cytometry (suitable as primary antibody in staining for FACS 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.

UniProt ID

P05362

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)