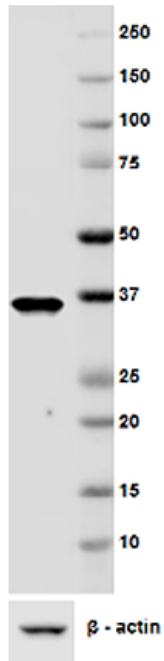


Lymphatic endothelial marker (human) monoclonal antibody (D2-40)

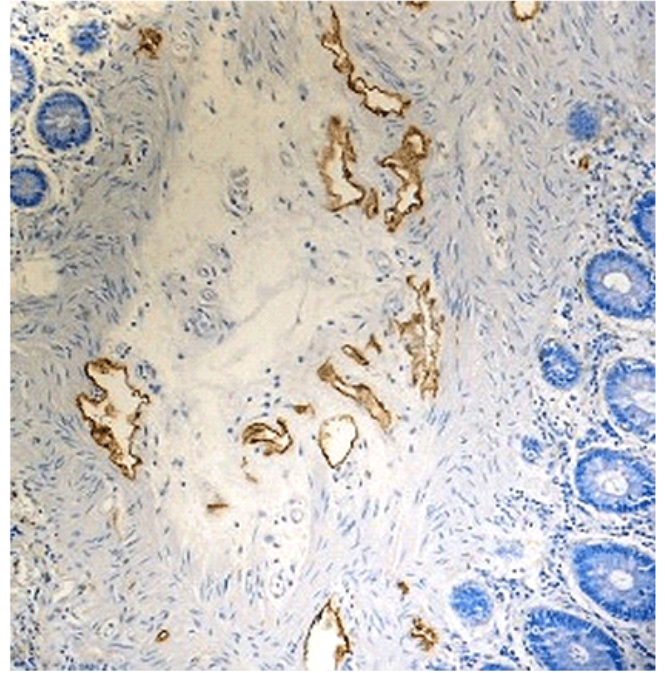
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of cell lysates from HeLa using Podoplanin Mouse primary antibody and HRP Goat anti-Mouse secondary antibody.



IHC staining of purified anti-Podoplanin (Lymphatic Endothelial Marker) antibody (clone D2-40) on formalin-fixed paraffin embedded human colon tissue.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Podoplanin, Aggrus, Glycoprotein 36, T1A

Application IHC, WB

Clone D2-40

Formulation Liquid. In Phosphate-buffered solution, pH 7.2.

Host Mouse

Immunogen This antibody was developed using M2A.

Isotype IgG1

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions Each lot of this antibody is quality control tested by immunohistochemistry. For immunohistochemistry on formalin-fixed paraffin-embedded tissue, a concentration range of 0.5 – 5.0 µg/ml is suggested. For Western blotting, the suggested use of this reagent is 2.0 – 5.0 µg per ml. It is recommended that the reagent be titrated for optimal performance for each application.

Species Reactivity Human

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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