

Vinculin (human) monoclonal antibody (FB11)

Vinculin is associated with the cytoplasmic aspect of adhesion points and acts as a link between the ends of actin microfilaments and the plasma membrane. Vinculin may also be involved in the transmembrane induction of actin-bundle formation.

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

Ordering Information [Order Online »](#)

BML-VG6110-0100	100µl
-----------------	-------

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	IHC, WB
Clone	FB11
Formulation	Liquid. In PBS containing 1.0% BSA and 0.09% sodium azide.
Host	Mouse
Immunogen	Purified vinculin from human smooth muscle cells.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Species Reactivity	Human
UniProt ID	P18206
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-be@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