

Fibronectin (ED-B domain) Recombinant monoclonal antibody (A2) (Human IgG1κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS633-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Extra domain-B containing fibronectin
Application	ELISA, IHC (FS), WB
Clone	A2
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	2D-PAGE gel isolated human Extra-Domain containing fibronectin from lysate of human melanoma COLO-38 cells.
Isotype	IgG1k
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Technical Info / Product Notes	This full-length, reformatted human antibody was made using the variable domain sequences of the original Human scFv format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID P02751



Worry-free Guarantee

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +32 3 466 0420
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
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

This antibody is covered by our Worry-Free Guarantee