

Fibroblast activation protein (human) monoclonal antibody (F11-24)

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

Ordering Information

[Order Online »](#)

ALX-804-327-C100	100µg
------------------	-------

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	FAP
Application	Flow Cytometry, IHC (FS), IP
Application Notes	Not suitable for Immunohistochemistry (paraffin sections).
Clone	F11-24
Crossreactivity	Cross-reacts with a subset of endocrine cells in the pancreatic islets, presumably glucagon-producing A cells.
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Human FAP (fibroblast activation protein).
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (suitable for stainings on unfixed, acetone-fixed or ethanol-fixed frozen sections; acetone-fixed frozen sections showed optimal staining results)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human

Technical Info / Product Notes	FAP is a cell surface glycoprotein of reactive human fibroblasts. Among the normal adult organs tested, MAb to FAP does not show immunoreactivity in immunohistological studies with either resting fibroblasts (fibrocytes) of the normal organ mesenchyme or the parenchymal cells. The single exception noted is a subset of endocrine cells in the pancreatic islets, presumably glucagon-producing A cells.
UniProt ID	Q12884
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)