

# FABP4 recombinant monoclonal antibody (001)

Fatty Acid Binding Protein 4 (FABP4) is primarily expressed in adipocytes and macrophages, playing a significant role in lipid metabolism, inflammation, and insulin sensitivity. It is widely studied in metabolic disorders, obesity, diabetes, and cardiovascular diseases as a potential biomarker and therapeutic target.

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

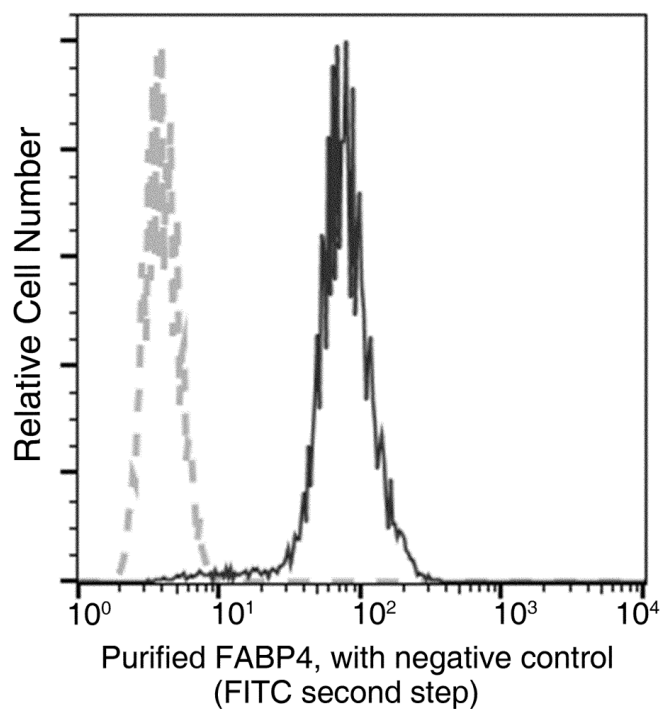
## Ordering Information

[Order Online »](#)

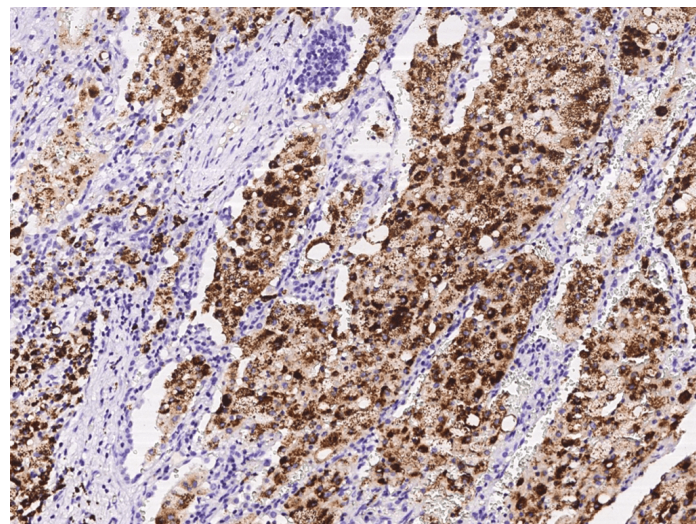
ENZ-ABS1032-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human FABP4 expression on Hela cells. The cells were treated according to manufacturer's manual (BD Pharmingen™ Cat. No. 554714), stained with purified anti-Human FABP4, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunochemical staining of human FABP4 in human lung cancer with rabbit monoclonal antibody (1:2000, formalin-fixed paraffin embedded sections).

# Handling & Storage

Handling	Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Fatty acid binding protein 4, Adipocyte fatty acid binding protein, ALBP, Adipocyte lipid binding protein
Application	Flow Cytometry, IHC (PS)
Clone	001
Crossreactivity	No cross-reactivity in ELISA with human FABP6.
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free. Carrier-free.
Host	Rabbit
Immunogen	Recombinant human FABP4 / ALBP / A-FABP.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:1'000-1:4'000)Flow Cytometry (1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)