

# Complement component C5a polyclonal antibody

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

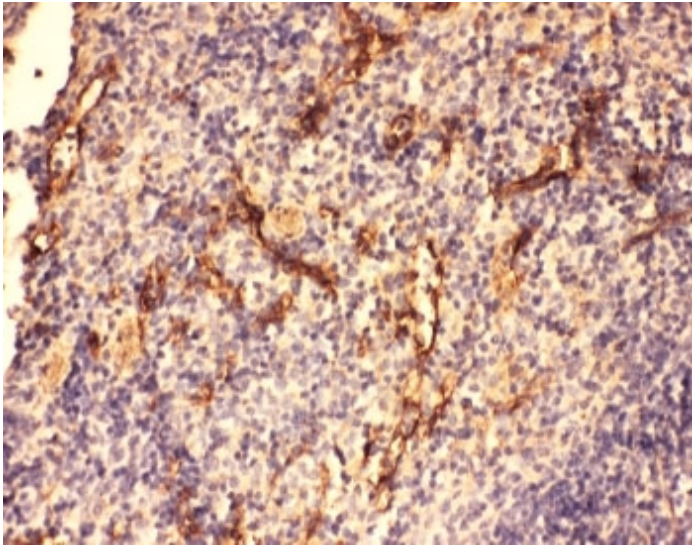
## Ordering Information

[Order Online »](#)

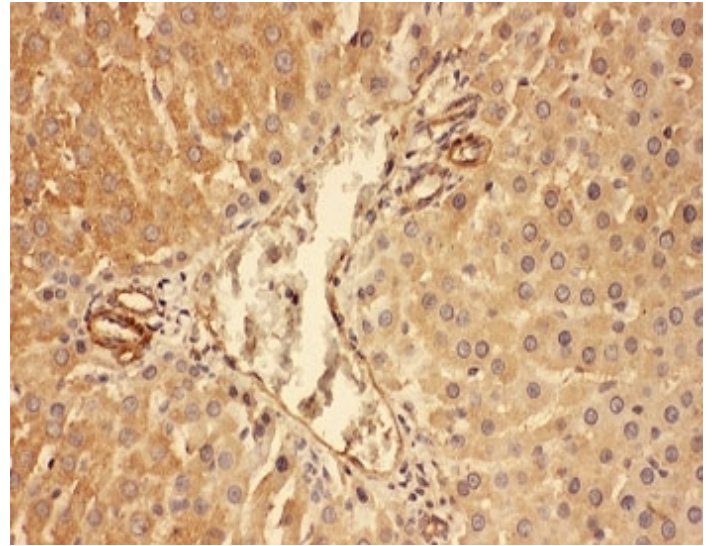
ENZ-ABS874-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

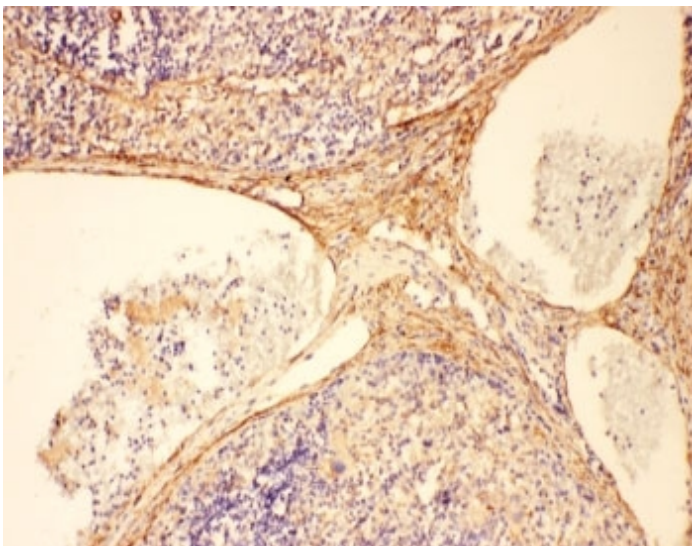
[View Online »](#)



IHC-P: C5a antibody testing of mouse spleen tissue



IHC-P: C5a antibody testing of rat liver tissue



IHC-P: C5a antibody testing of rat spleen tissue



Western blot testing of C5a antibody and Lane 1: rat liver; 2: mouse liver; 3: (m) NIH3T3; 4: (m) HEPA; 5: (r) PC12 lysate; Predicted/Observed size: 115KD

## Handling & Storage

<b>Use/Stability</b>	After reconstitution, the C5a antibody can be stored for up to one month at +4°C. For long-term, aliquot and store at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Application</b>	IHC (PS), WB
<b>Formulation</b>	Lyophilized from 1X PBS containing 2.5% BSA and 0.025% sodium azide/thimerosal.
<b>Host</b>	Rabbit
<b>Immunogen</b>	Amino acids DLQLLHQKVVEEQAAYKH were used as the immunogen for this C5a antibody.
<b>Isotype</b>	IgG
<b>Purity Detail</b>	Antigen affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Immunohistochemistry (FFPE) (0.5-1µg/ml)Western blot (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Mouse, Rat
<b>Worry-free Guarantee</b>	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-  
fr@enzolifesciences.com](mailto:info-fr@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)