

# Complement component C3a/C3a (desArg) monoclonal antibody (K13/16)

Complement C3a is an anaphylatoxin of 77 amino acid residues released by the action of the C3 convertases on the N-terminal of the alpha chain of C3. It is rapidly inactivated by serum carboxypeptidase N which removes the C-terminal arginine residue generating C3a (desArg).

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

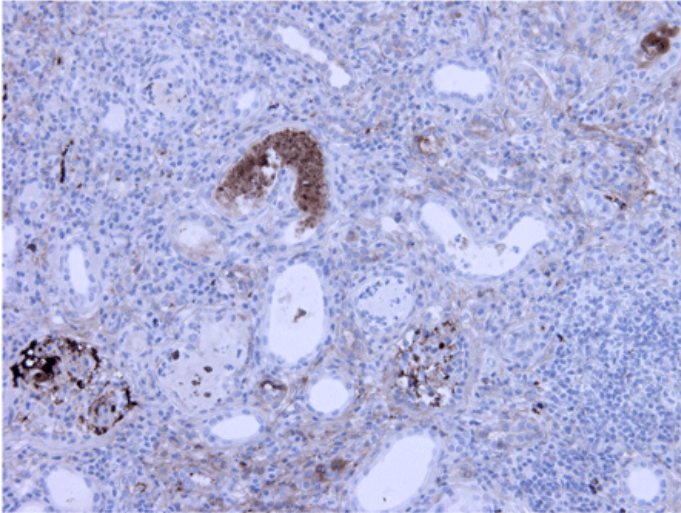
## Ordering Information

[Order Online »](#)

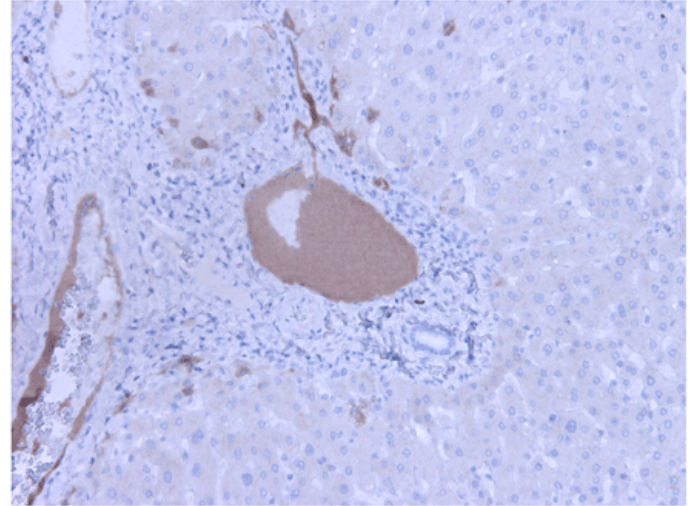
BPD-GAU-013-16-04	400µg
-------------------	-------

## Manuals, SDS & CofA

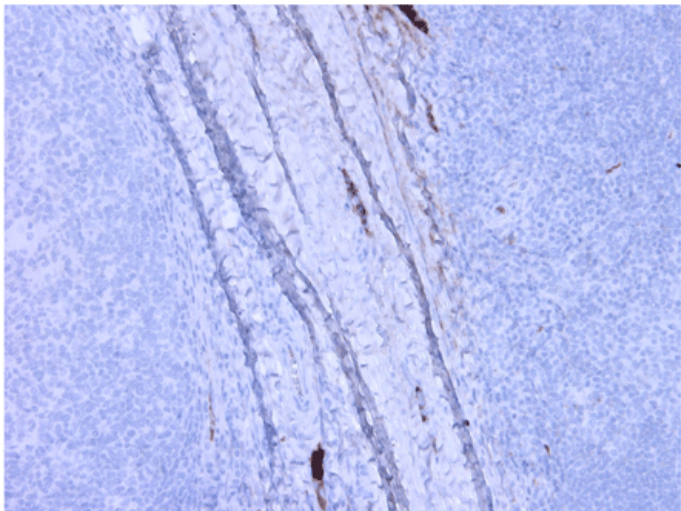
[View Online »](#)



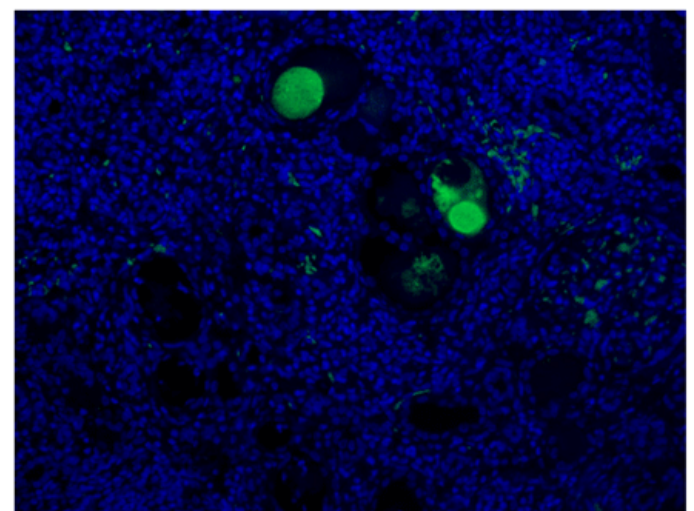
IHC staining of human kidney with BPD-GAU-013-16.  
Strong staining of veins of nephritis kidney was seen.  
Dilution: 1:150.



IHC staining of human liver with BPD-GAU-013-16.  
Strong staining of plasma of veins in liver tissue.  
Dilution: 1:150.



IHC staining of human tonsil with BPD-GAU-013-16.  
Plasma in lumen of veins is stained strongly in tonsil.  
Dilution: 1:150.



IF staining in human kidney with BPD-GAU-013-16.  
Strong staining of plasma of nephritis kidney was seen.  
Dilution: 1:75.

## Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	ELISA, IF, IHC, WB
Clone	K13/16
Crossreactivity	Does not cross-react with C4a or C5a.
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Human complement component C3a.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Human
Specificity	Recognizes C3, C3a and C3a (desArg).
UniProt ID	P01024
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)