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.

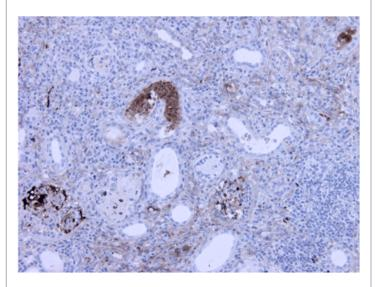
Ordering Information

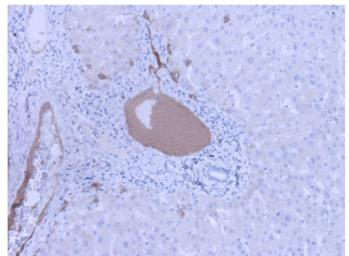
Order Online »

| BPD-GAU-013-16-1 | 1mg |
|-------------------|-------|
| BPD-GAU-013-16-04 | 400µg |

Manuals, SDS & CofA

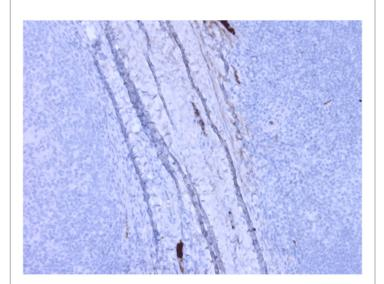
View Online »

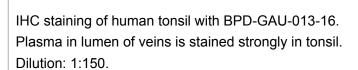


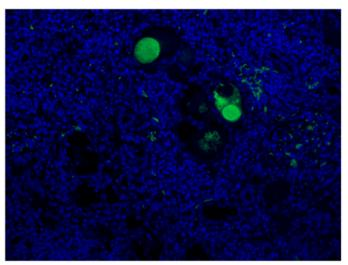


IHC staining of human kidney with BPD-GAU-013-16. Strong staining of veins of nephritis kidney was seen. Dilution: 1:150.

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







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.

lgG1κ

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 Worry-Free Guarantee

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 info-

usa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com