Complement component C5a polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

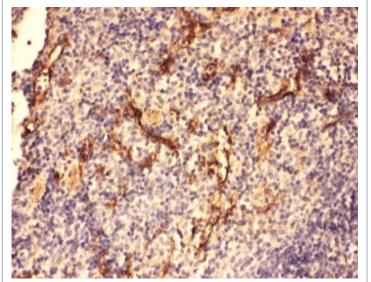
Order Online »

ENZ-ABS874-0100

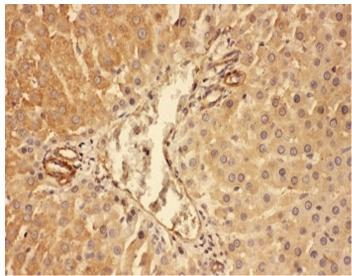
100µg

Manuals, SDS & CofA

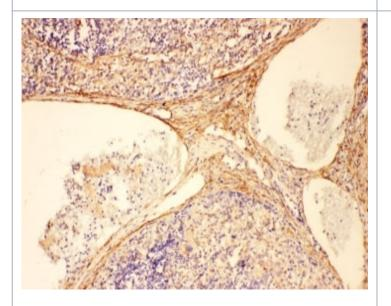
View Online »



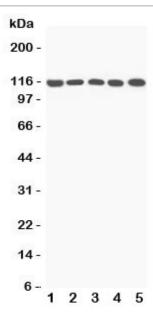
IHC-P: C5a antibody testing of mouse spleen tissue



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

Use/Stability 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.

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Formulation Lyophilized from 1X PBS containing 2.5% BSA and

0.025% sodium azide/thimerosal.

Host Rabbit

Immunogen Amino acids DLQLLHQKVEEQAAKYKH were used as the

immunogen for this C5a antibody.

Isotype IgG

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions 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.

Species Reactivity Mouse, Rat

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•



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