Complement component C5 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

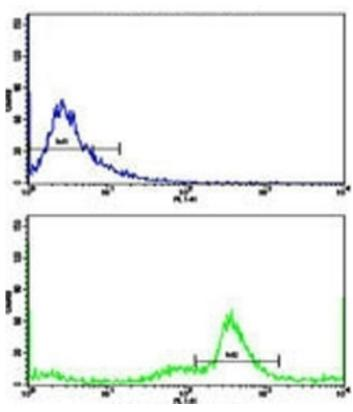
Order Online »

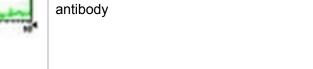
ENZ-ABS873-0400

0.4ml

Manuals, SDS & CofA

View Online »





IHC analysis of FFPE human breast carcinoma with C5

Flow cytometric analysis of MDA-231 cells using C5 antibody (green) compared to a negative control (blue). FITC-conjugated goat-anti-rabbit secondary Ab was used for the analysis.

Western blot analysis of C5 antibody and HL-60 lysate

Handling & Storage

Handling Aliquot the C5 antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, Flow Cytometry, IHC, WB

Formulation Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.

Host Rabbit

Immunogen Human C5 (aa 253-281).

Isotype Ig

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin

sections) (1:50-1:100)Western blot (1:1,000)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P01031

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

.

Last modified: July 12, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@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