LYPD3 recombinant monoclonal antibody (213)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

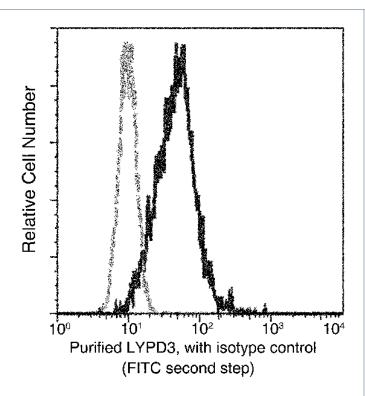
Order Online »

ENZ-ABS983-0100

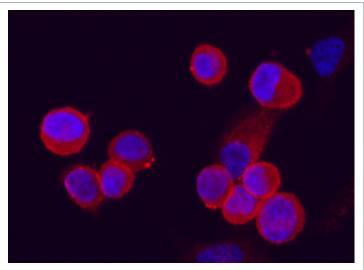
100µl

Manuals, SDS & CofA

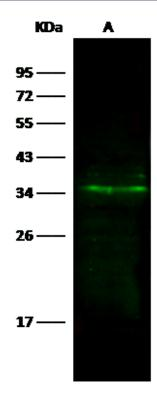
View Online »



Flow cytometric analysis of Human LYPD3 expression on PC-3 cells. Cells were stained with purified anti-Human LYPD3, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of Human LYPD3 in PC3 cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-Human LYPD3 monoclonal antibody (1:60) at 37 1 hour. Then cells were stained with the Alexa Fluor® 594-conjugated goat Anti-rabbit IgG secondary antibody (red) and counterstained with DAPI (blue). Positive staining was localized to cells membrane.



Anti-LYPD3 rabbit monoclonal antibody at 1:500 dilution. Lane A: Hela Whole Cell Lysate:
Lysates/proteins at 30 µg per lane. Secondary Goat
Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution.
Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:36 kDa.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name C4.4A, Ly6/PLAUR domain-containing protein 3

Application ELISA, Flow Cytometry, ICC, IF, WB

Clone 213

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human C4.4A / LYPD3 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Immunocytochemistry (1:20-

1:100)Immunofluorescence (1:20-1:100)Flow Cytometry (1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

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

September 77, 20 Suppean Sales Office ENZO LIFE SCIENCES
Phone: 800.942.0430 (ELS) AG
infousa@enzolifesciences.com
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