

LYPD3 recombinant monoclonal antibody (213)

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

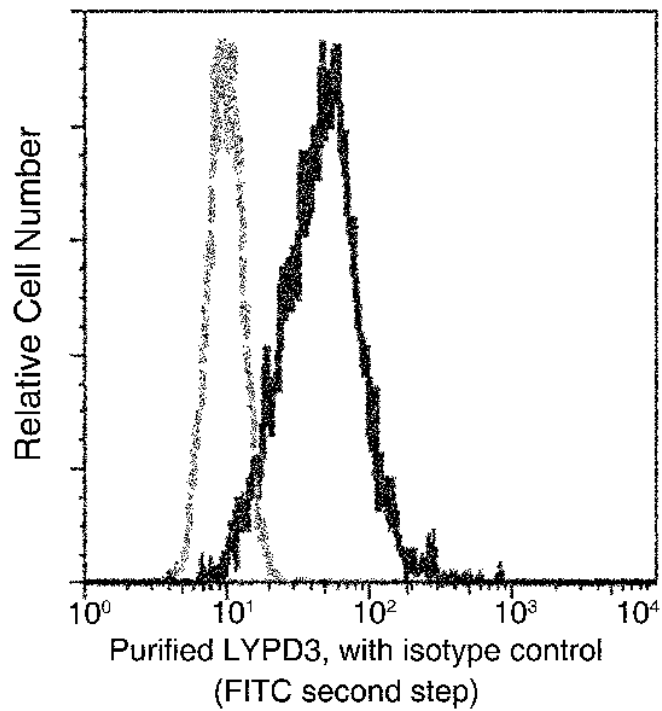
Ordering Information

[Order Online »](#)

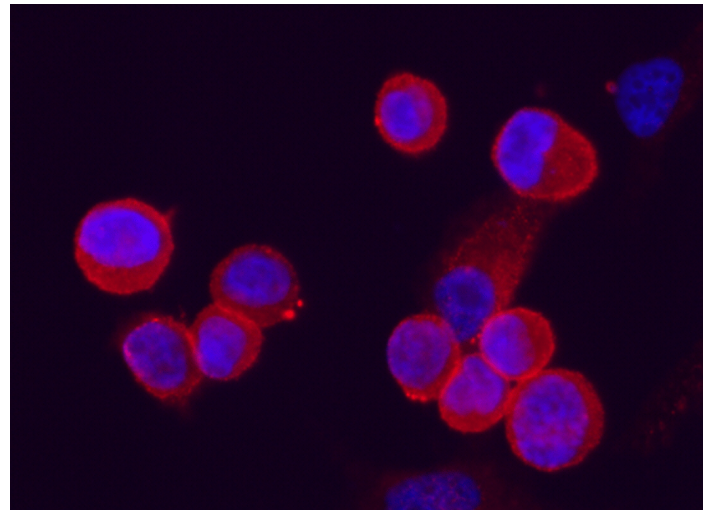
ENZ-ABS983-0100	100µl
-----------------	-------

Manuals, SDS & CofA

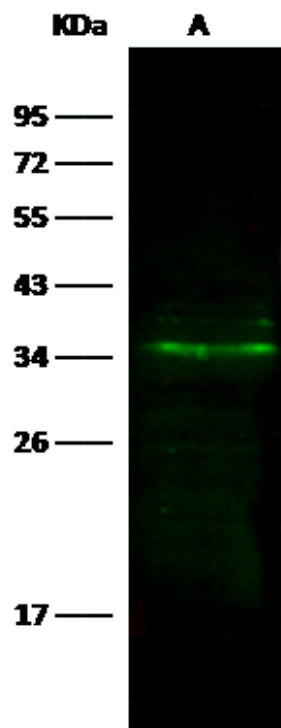
[View Online »](#)



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 °C for 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 μ g 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](#).



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-
fr@enzolifesciences.com](mailto:info-fr@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)