

OLFM4 monoclonal antibody (12)

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

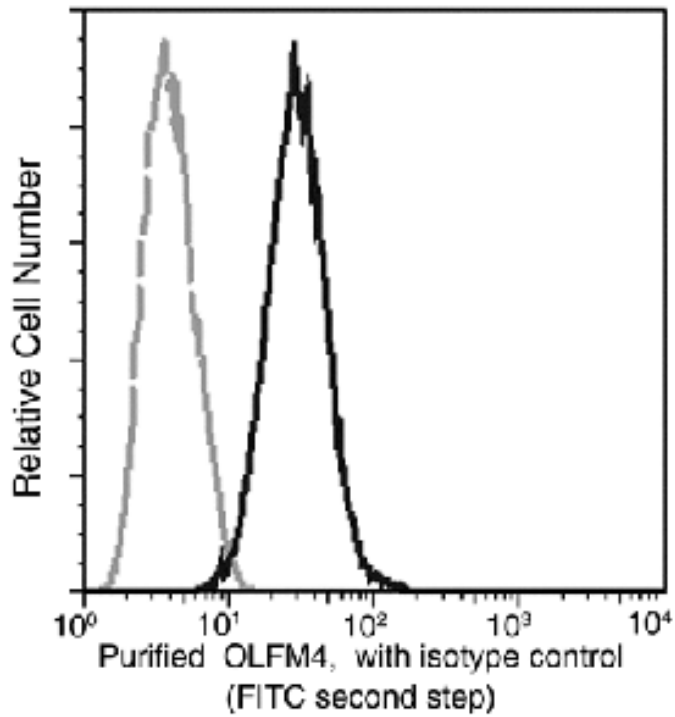
Ordering Information

[Order Online »](#)

ENZ-ABS981-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human OLFM4 expression in DU145 cells. The cells were treated and stained with Purified Mouse anti-OLFM4 (11639-MM12), 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.

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 Olfactomedin-4, OLM4, G-CSF-stimulated clone 1 protein, hGC-1, GW112

Application ELISA, Flow Cytometry

Clone 12

Formulation Liquid. 0.2 µm filtered solution in PBS with 10% Trehalose, pH 7.0. Preservative free.

Host Mouse

Immunogen Recombinant human OLFM4 / GW112 protein.

Isotype IgG1

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000)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: 800 642 0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EU) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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