

Langerin (human) monoclonal antibody (DCGM4/122D5) (Alexa Fluor[®] 488)

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

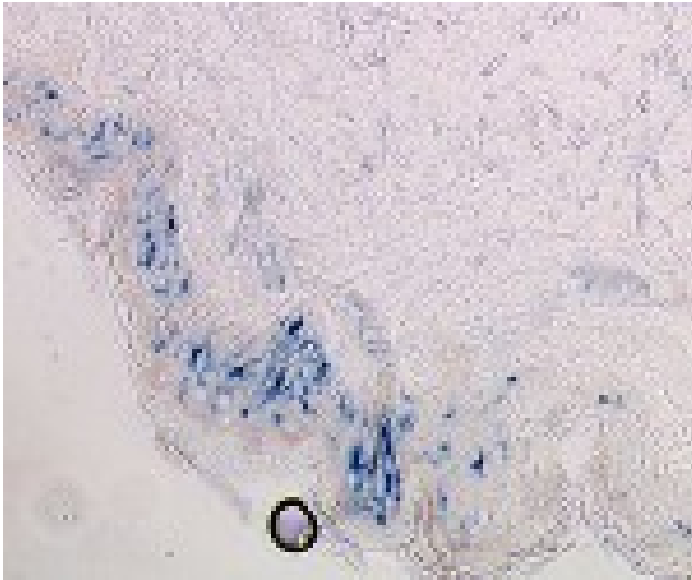
Ordering Information

[Order Online »](#)

ENZ-ABS238-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human skin frozen section with Langerin (human) mAb, clone DCGM4/122D5.

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD207

Application Flow Cytometry, IF, IHC (FS)

Clone DCGM4/122D5

Formulation Liquid. In PBS containing 50% glycerol.

Host Mouse

Immunogen In vitro-derived human DCs from CD34+ cells in GM-CSF+TNF- α .

Isotype IgG1

Purity Detail Purified using QMA Hyper D ion exchange chromatography

Species Reactivity Dog, Human

Specificity Recognizes an extracellular epitope in carbohydrate recognition domain of Langerin.

UniProt ID Q9UJ71

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