

Langerin (human) monoclonal antibody (DCGM4/122D5)

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

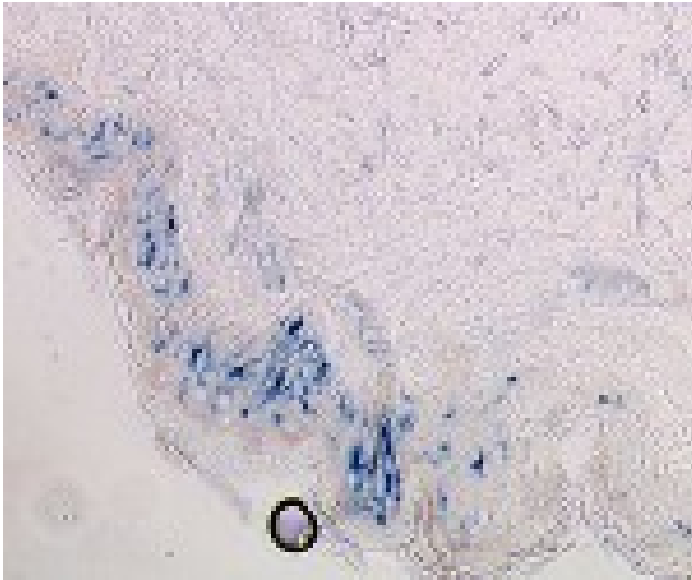
Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS242-0100 | 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, IHC (FS), Internalization |
| Clone | DCGM4/122D5 |
| Formulation | Liquid. In TRIS-NaCl, pH 8. |
| 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](#)

.



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@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