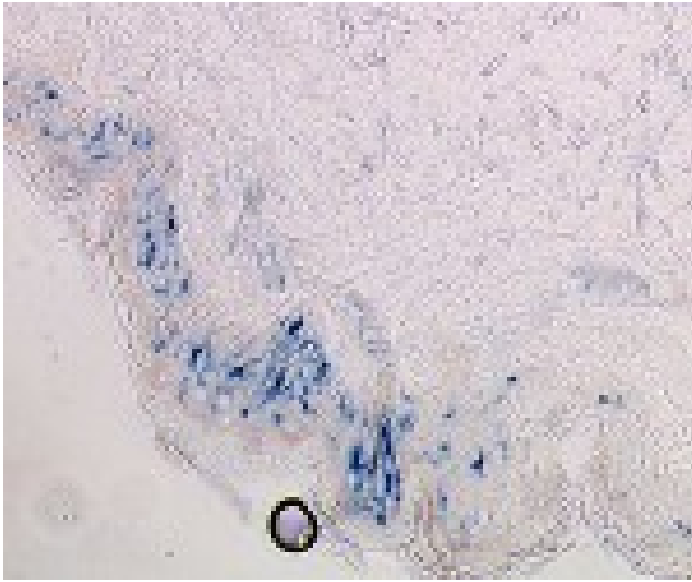


# Langerin (human) monoclonal antibody (DCGM4/122D5) (Alexa Fluor<sup>®</sup> 546)

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

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	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- $\alpha$ .
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  
Last modified: **May 29, 2024**  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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