

# IL-3 receptor $\alpha$ (human) monoclonal antibody (107D2.08)

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

Manuals, SDS & CofA

[View Online »](#)

## 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** CD123, IL-3RA, Interleukin-3 receptor  $\alpha$

**Application** Flow Cytometry, IHC (FS), IHC (PS), IP

**Clone** 107D2.08

**Formulation** Liquid. In TRIS-NaCl, pH 8.

**Host** Mouse

**Immunogen** Sorted human tonsillar pDC.

**Isotype** IgG1

**Purity Detail** Purified using QMA Hyper D ion exchange chromatography

**Species Reactivity** Dog, Human

**Specificity** Recognies an epitope in the extracellular domain of CD123.

**UniProt ID** P26951

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



ENZO LIFE SCIENCES,  
**Last modified May 29, 2024**  
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)