

IL-13 receptor (human) monoclonal antibody (GM-1C8)

IL-13 is an immunoregulatory cytokine secreted predominantly by activated TH2 cells, as well as mast cells and basophils, and is a key mediator in the pathogenesis of allergic inflammation. The human IL-13 receptor (IL-13R) is a heterodimer composed of the interleukin-4 receptor α chain (IL-4R α) and an IL-13 binding protein, IL-13R α 1. A second IL-13 binding protein, IL-13R α 2, has also been identified, which is believed to act as a "decoy" receptor. IL-13 receptors are expressed on human B cells, basophils, eosinophils, mast cells, endothelial cells, fibroblasts, monocytes, macrophages, respiratory epithelial cells, and smooth muscle cells. However, functional IL-13 receptors have not been demonstrated on human or mouse T cells.

As with other cytokine receptors, the cytoplasmic tails of IL-4 and IL-13 receptor subunits associate with tyrosine kinases of the Janus family (JAK 1

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-506-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

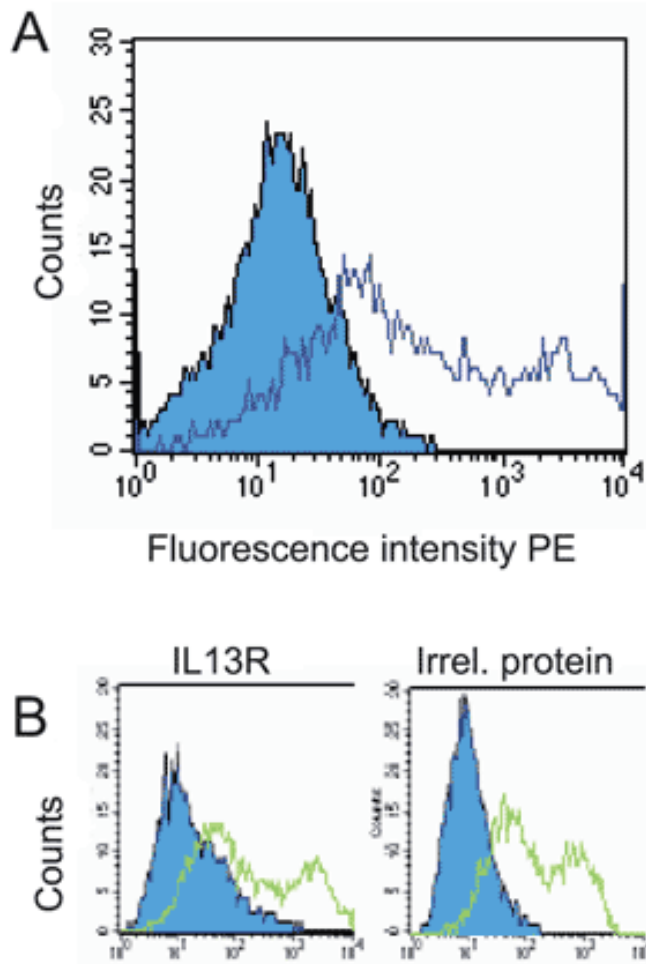


Figure: Specificity testing of GM-1C8. BOSC cells were transiently transfected with an expression vector for the extracellular domain of IL-13Ra1 as well as an irrelevant protein. Expression of the constructs was tested with an anti-tag monoclonal antibody (B, open curves), an irrelevant monoclonal antibody served as negative control (filled curves). For specificity testing (A), GM-1C8 hybridoma supernatant was tested on both the IL-13Ra1 and the irrelevant protein transfectants. A positive signal was obtained only with IL-13Ra1-transfected cells.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	IL-13R, CD213, Interleukin-13 receptor
Application	Flow Cytometry
Clone	GM-1C8
Formulation	Liquid. Antibody in PBS, pH 7.2, containing 0.01% sodium azide.
Host	Mouse
Immunogen	Vector containing the cDNA of the extracellular domain of human IL-13Rα1 (interleukin-13 receptor α1 chain).
Isotype	IgG1
Quality Control	Routinely tested by flow cytometry on BOSC cells transiently transfected with a IL-13Rα1 expression vector.
Source	Purified by affinity chromatography from cell culture supernatant.
Species Reactivity	Human
UniProt ID	P78552 (subunit α1), Q14627 (subunit α2)
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)