

LT β R (mouse) monoclonal antibody (5G11)

LT β R is involved in peripheral lymphoid tissue organogenesis and function. The LT α /LT β R receptor system may also have some function in early embryogenesis. LT β R is a type I transmembrane glycoprotein and member of the TNF receptor superfamily. It is a 61kDa receptor for the mouse LT α 1 β 2 and LIGHT. Northern blot analysis of tissues from adult mice showed that expression levels of LT β R mRNA were strong in lung, liver, and kidney, moderate in heart and testes, but weak in brain, thymus, spleen, and lymph nodes. It was reported that LT β R is expressed on stromal cells, monocytes, and certain embryonic epithelial layers. LT β R is involved in peripheral lymphoid tissue organogenesis and function. The LT α /LT β R receptor system may also have some function in early embryogenesis.

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

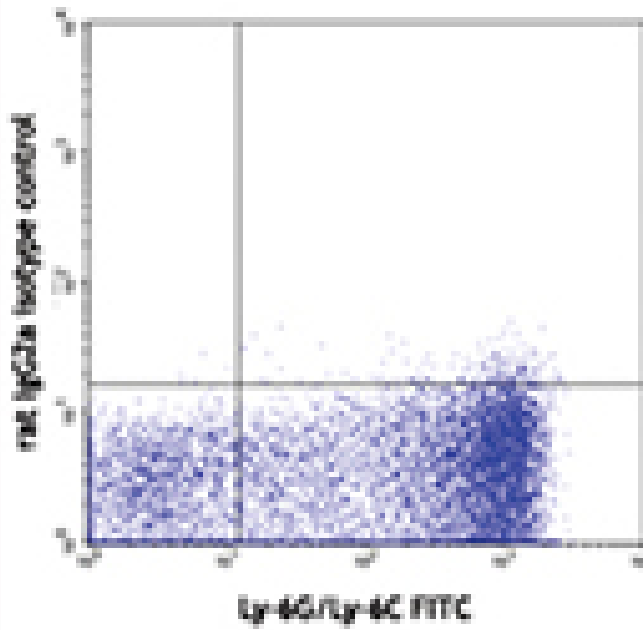
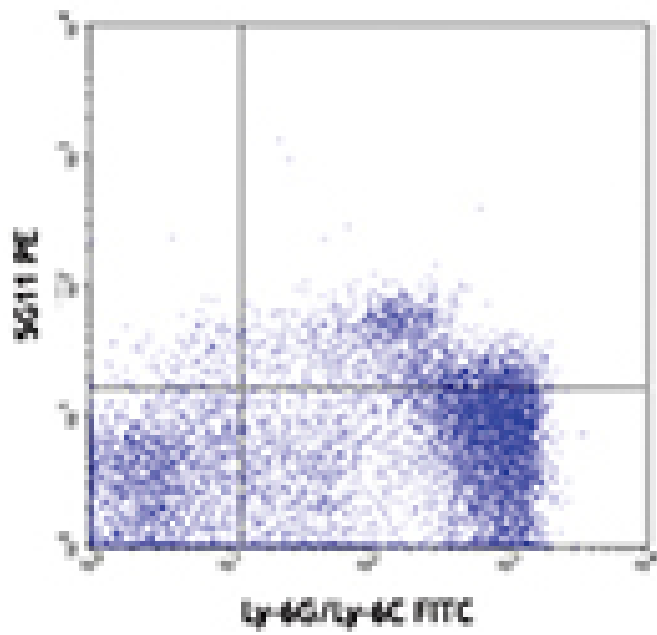
Ordering Information

[Order Online »](#)

ALX-804-903-0100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of C57BL/6 bone marrow cells stained with purified LT β R (mouse), mAb (5G11) (Prod. No. ALX-804-903) conjugated to PE (Upper panel) or rat IgG2a isotype control PE (Lower panel) and Ly-6G/Ly-6C FITC.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name

Lymphotoxin- β receptor, TNFRSF 3, Tumor necrosis factor receptor 2 related protein, Tumor necrosis factor C receptor, Tumor necrosis factor receptor superfamily member 3 [precursor]

Application

Flow Cytometry

Clone

5G11

Formulation

Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Gene/Protein Identifier

17000 (Entrez GeneID)

Host

Rat

Immunogen

Extracellular domain of mouse LT β R (lymphotoxin β receptor) expressed in E. coli.

Isotype

IgG2a

Purity Detail

Affinity purified.

Species Reactivity

Mouse

UniProt ID

P50284

Worry-free Guarantee

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

.