

# LT $\beta$ R (mouse) monoclonal antibody (5G11)

LT $\beta$ R is involved in peripheral lymphoid tissue organogenesis and function. The LT $\alpha$ /LT $\beta$ R receptor system may also have some function in early embryogenesis. LT $\beta$ R is a type I transmembrane glycoprotein and member of the TNF receptor superfamily. It is a 61kDa receptor for the mouse LT $\alpha$ 1 $\beta$ 2 and LIGHT. Northern blot analysis of tissues from adult mice showed that expression levels of LT $\beta$ 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 $\beta$ R is expressed on stromal cells, monocytes, and certain embryonic epithelial layers. LT $\beta$ R is involved in peripheral lymphoid tissue organogenesis and function. The LT $\alpha$ /LT $\beta$ R receptor system may also have some function in early embryogenesis.

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

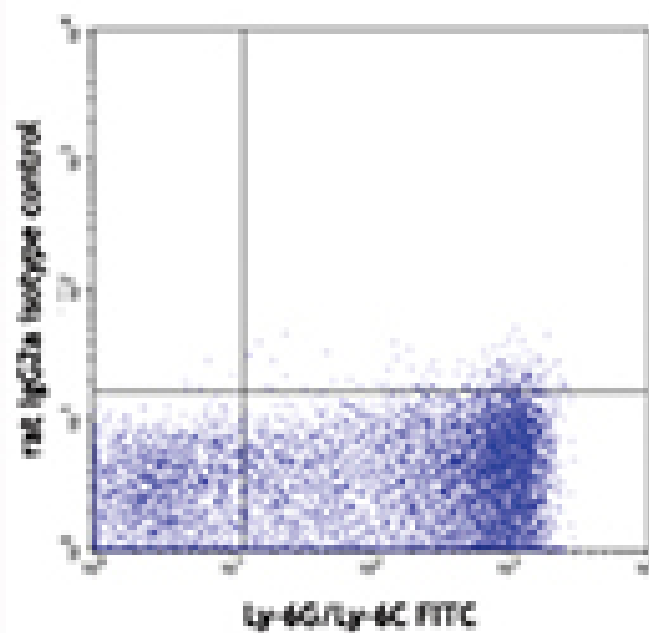
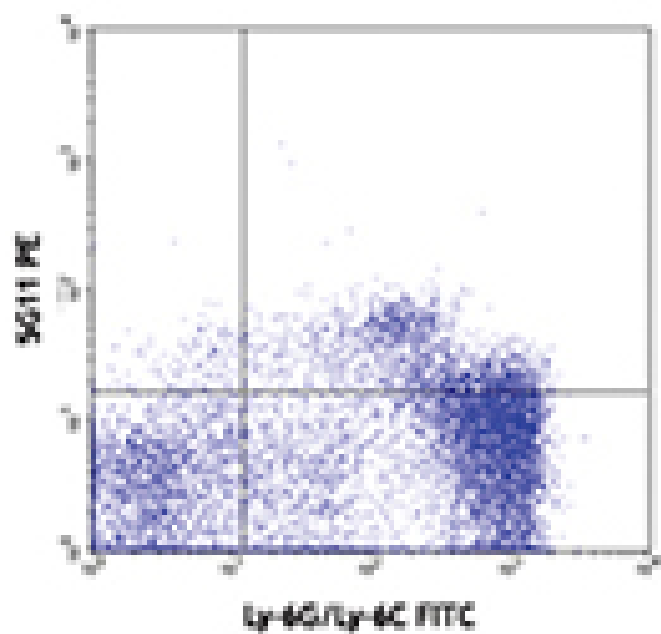
## Ordering Information

[Order Online »](#)

ALX-804-903-0100	100 $\mu$ g
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of C57BL/6 bone marrow cells stained with purified LT $\beta$ 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- $\beta$  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 $\beta$ R (lymphotoxin  $\beta$  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](#).



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)