

TLR7 monoclonal antibody (4F4)

TLR7 is one of toll-like receptors which are type I transmembrane proteins involved in innate immunity by recognizing microbial conserved structures. These proteins are characterized by multiple copies of leucine rich repeats in the extracellular domain and IL-1 receptor motif in the cytoplasmic domain, and act through adaptor molecules such as MyD88 and TIRAP to activate various kinases and transcription factors.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	toll-like receptor 7
Application	ELISA, Flow Cytometry, ICC, IF
Clone	4F4
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
Host	Mouse
Immunogen	Recombinant human TLR7 (aa 451-500) purified from E. coli.
Isotype	IgG1κ
Purity Detail	Protein G affinity purified
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	Q9NYK1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee