

Tlr7 Polyclonal Antibody

(Catalog # A53663)

Background

Key component of innate and adaptive immunity. TLRs (Toll-like receptors) control host immune response against pathogens through recognition of molecular patterns specific to microorganisms. TLR7 is a nucleotide-sensing TLR which is activated by single-stranded RNA. Acts via MYD88 and TRAF6, leading to NF-kappa-B activation, cytokine secretion and the inflammatory response

Description

Tlr7 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P58681

Purification

Protein G purified

Immunogen

Recombinant Mouse Toll-like receptor 7 protein (27-348AA)

Storage

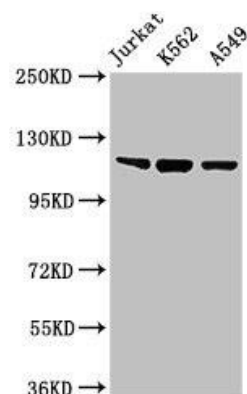
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Tlr7

Application

ELISA, WB; Recommended dilution: WB:1:500-1:2000



Western Blot

Positive WB detected in: Jurkat whole cell lysate, K562 whole cell lysate, A549 whole cell lysate

All lanes: Tlr7 Polyclonal Antibody at 2ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 122 KDa

Observed band size: 122 KDa