

LGP2 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat LGP2 (CT)

Instruction Manual

Catalog Number	PK-AB718-4351
Synonyms	LGP2 Antibody: DHX58, likely ortholog of mouse D11lgp2
Description	Anti-viral innate immune responses are triggered by pathogen-associated molecular patterns (such as the accumulation of intracellular nucleic acids resulting from virus infections) and represent the first line of defense against numerous infectious organisms. The Toll-like receptors (TLR) 3, 7, 8, and 9 are expressed in immune cells and function as transmembrane pattern recognition receptors to detect foreign nucleic acids. Other proteins that play similar roles, such as RIG-1 and MDA5 are members of a CARD-helicase family and are expressed in the cytoplasm. A third protein, LGP2, is similar to RIG-1 and MDA5, except for lacking a homologous CARD domain. It is thought to act as an element of negative-feedback regulation of intracellular antiviral signaling. At least four isoforms of LGP2 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	LGP2 antibody was raised in rabbits against a 12 amino acid peptide from near the carboxy terminus of human LGP2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. LGP2 antibody can be used for detection of LGP2 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Akira S, Uematsu S, and Takeuchi O. Pathogen recognition and innate immunity. <i>Cell</i> 2006; 124:783-801. Iwasaki A and Medzhitov R. Toll-like receptor control of the adaptive immune responses. <i>Nat. Immunol.</i> 2004; 5:987-95. Yoneyama M, Kikuchi K, Matsumoto K, et al. The RNA helicase RIG-I has an essential function in double-stranded RNA-induced innate antiviral responses. <i>Nat. Immunol.</i> 2004; 5:730-7. Andrejeva J, Childs KS, Young DF, et al. The V proteins of paramyxoviruses bind the IFN-inducible RNA helicase, mda-5, and inhibits its activation of the IFN-beta promoter. <i>Proc. Natl. Acad. Sci. USA</i> , 2004; 101:17264-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4351P; LGP2 Peptide Cat.No. PK-AB718-1465; Rat Kidney Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.