

Cathelicidin antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Cathelicidin (CAP18, FALL39, LL37)

Instruction Manual

Catalog Number	PK-AB718-4429
Synonyms	Cathelicidin Antibody: Cathelicidin, Cathelicidin antimicrobial peptide, Cationic antimicrobial protein, CAP18, FALL39, hCAP-18, HSD26, LL37
Description	One component of host defense at mucosal surfaces is epithelial-derived antimicrobial peptides. Cathelicidins are one family of antimicrobial peptides characterized by conserved pro-peptide sequences that have been identified in epithelial tissues and some myeloid cells of humans and animals. LL-37/hCAP-18 is the only Cathelicidin found in humans and is expressed in inflammatory and epithelial cells. The presence of these molecules is essential for defense against invasive bacterial infection in skin. Besides their direct antimicrobial function, Cathelicidins have multiple roles in mediating innate and adaptive immunity, such as endotoxin neutralizing, angiogenesis, wound healing and promoting neutrophil chemotaxis and mast cell recruitment. Finally, Cathelicidin antimicrobial peptides qualify as prototypes of innovative drugs that may be used to treat infection and/or modulate the immune response.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Cathelicidin antibody was raised in rabbits against a 17 amino acid peptide near the carboxy terminus of the human Cathelicidin.
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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Cathelicidin antibody can be used for detection of Cathelicidin by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Zaiou M and Gallo RL. Cathelicidins, essential gene-encoded mammalian antibiotics. <i>J. Mol. Med.</i> 2002; 80:549-61. Agerberth B, Gunne H, Odeberg J, et al. FALL-39, a Putative Human Peptide Antibiotic, is Cysteine-Free and Expressed in Bone Marrow and Testis. <i>Proc. Natl. Acad. Sci. USA</i> 1995; 92:195-9. Nizet V, Ohtake T, Lauthe X, et al. Innate antimicrobial peptide protects the skin from invasive bacterial infection. <i>Nature</i> 2001; 414:454-7. Koczulla R, von Degenfeld G, Kupatt C, et al. An angiogenic role for the human peptide antibiotic LL-37/hCAP-18. <i>J. Clin. Invest.</i> 2003; 111:1665-72.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4429P; Cathelicidin Peptide Cat.No. PK-AB718-1306; Human Spleen Tissue Lysate

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