

Instruction Manual

Catalog Number	PK-AB718-3665P
Quantity	50 µg
Source	17 amino acids near the carboxy terminus of human BAP31
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store BAP31 peptide at -20°C, stable for one year.
Application	BAP31 peptide is used for blocking the activity of BAP31 antibody.
References	<p>Kim KM, Adachi T, Nielsen PJ, et al. Two new proteins preferentially associated with membrane immunoglobulin D. <i>EMBO J.</i> 1994; 13:3793-800.</p> <p>McGettrick AF and O'Neill LAJ. The expanding family of MyD88-like adaptors in Toll-like receptor signal transduction. <i>Mol. Imm.</i> 2004; 41:577-82.</p> <p>Takeuchi O, Kawai T, Sanjo H, et al. TLR6: A novel member of an expanding Toll-like receptor family. <i>Gene</i> 1999; 231:59-65. 5. Ozinsky A, Underhill DM, Fontenot JD, et al. The repertoire for pattern recognition of pathogens by the innate immune system is defined by cooperation between toll-like receptors. <i>Proc. Natl. Acad. Sci. USA</i> 2000; 97:13766-71. 6. Into T, Kiura K, Yasuda M, et al. Stimulation of human Toll-like receptor (TLR) 2 and TLR6 with membrane lipoproteins of <i>Mycoplasma fermentans</i> induces apoptotic cell death after NF-kappa B activation. <i>Cell Microbiol.</i> 2004; 6:187-99.</p> <p>Shen J, Chen X, Hendershot L, et al. ER stress regulation of ATF6 localization by dissociation of BiP/GRP78 binding and unmasking of Golgi localization signals. <i>Dev. Cell</i> 2002; 3:99-111.</p> <p>Ozinsky A, Underhill DM, Fontenot JD, et al. The repertoire for pattern recognition of pathogens by the innate immune system is defined by cooperation between toll-like receptors. <i>Proc. Natl. Acad. Sci. USA</i> 2000; 97:13766-71. Into T, Kiura K, Yasuda M, et al. Stimulation of human Toll-like receptor (TLR) 2 and TLR6 with membrane lipoproteins of <i>Mycoplasma fermentans</i> induces apoptotic cell death after NF-kappa B activation. <i>Cell Microbiol.</i> 2004; 6:187-99.</p>

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

09.10.2014

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

PromoCell GmbH
Sickingenstr. 63/65
69126 Heidelberg
Germany

North America
Deutschland
France
United Kingdom
Other Countries

1 – 866 – 251 – 2860 (toll free)
0800 – 776 66 23 (gebührenfrei)
0800 90 93 32 (ligne verte)
0800 – 96 03 33 (toll free)
+49 6221 – 649 34 0

Email: info@promokine.info
www.promokine.info

08/2014