

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3667P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino acids near the center of human BAP29
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store BAP29 peptide at -20°C, stable for one year.
<b>Application</b>	BAP29 peptide is used for blocking the activity of BAP29 antibody.
<b>References</b>	<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>Bertolotti A, Zhang Y, Hendershot LM, et al. Dynamic interaction of BiP and ER stress transducers in the unfolded-protein response. <i>Nat. Cell Biol.</i> 2000; 2:326-32.</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>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>Lee K, Tirasophon W, X Shen, et al. IRE1-mediated unconventional mRNA splicing and S2P-mediated ATF6 cleavage merge to regulate XBP1 in signaling the unfolded protein response. <i>Genes Dev.</i> 2002; 16:452-66.</p> <p>Miyoshi K, Katayama T, Imaizumi K, et al. Characterization of mouse Ire1 alpha: cloning, mRNA localization in the brain and functional analysis in a neural cell line. <i>Brain Res. Mol. Brain Res.</i> 2000; 85:68-76.</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](mailto:info@promokine.info)  
[www.promokine.info](http://www.promokine.info)

08/2014