

Instruction Manual

Catalog Number	PK-AB718-3029P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human Bmf
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 Bmf peptide at -20°C, stable for one year.
Application	Bmf peptide is used for blocking the activity of Bmf antibody.
References	<p>Puthalakath H, Villunger A, O'Reilly LA, Beaumont JG, Coultas L, Cheney RE, Huang DC, Strasser A. Bmf: a proapoptotic BH3-only protein regulated by interaction with the myosin V actin motor complex, activated by anoikis. <i>Science</i>. 2001;293(5536):1829-32</p> <p>Funke L, Dakoji S, Bredt DS. Membrane-Associated Guanylate Kinases Regulate Adhesion and Plasticity at Cell Junctions. <i>Annu. Rev. Biochem.</i> 2005; 74:219-45.</p> <p>Bilder D and Perrimon N. Localization of apical epithelial determinants by the basolateral PDZ protein Scribble. <i>Nature</i> 2000; 403:676-80.</p>

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