

Instruction Manual

Catalog Number	PK-AB718-2101P
Quantity	50 µg
Source	16 amino amino terminus of human GPR15
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 GPR15 peptide at -20°C, stable for one year.
Application	GPR15 peptide is used for blocking the activity of GPR15 antibody.
References	<p>Heiber M, Marchese A, Nguyen T, Heng HH, George SR, O'Dowd BF. A novel human gene encoding a G-protein-coupled receptor (GPR15) is located on chromosome 3. <i>Genomics</i> 1996;32:462-5.</p> <p>Liu X, Li P, Widlak P, Zou H, Luo X, Garrard WT, Wang X The 40-kDa subunit of DNA fragmentation factor induces DNA fragmentation and chromatin condensation during apoptosis. <i>Proc Natl Acad Sci USA</i> 1998;95:8461-6</p> <p>Inohara N, Koseki T, Chen S, Wu X, Nunez G. CIDE, a novel family of cell death activators with homology to the 45 kDa subunit of the DNA fragmentation factor. <i>EMBO J</i> 1998;17:2526-33</p>

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