

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-2051P |
| Quantity | 50 µg |
| Source | 17 amino acids near the carboxy terminus of human ADAM10 |
| 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 ADAM10 peptide at -20°C, stable for one year. |
| Application | ADAM10 peptide is used for blocking the activity of ADAM10 antibody. |
| References | <p>Rosendahl MS, Ko SC, Long DL, et al. Identification and characterization of a pro-tumor necrosis factor-alpha-processing enzyme from the ADAM family of zinc metalloproteases. <i>J Biol Chem</i> 1997;272:24588-93.</p> <p>Li P, Nijhawan D, Budihardjo I, Srinivasula SM, Ahmad M, Alnemri ES, Wang X. Cytochrome c and dATP-dependent formation of Apaf-1/caspase-9 complex initiates an apoptotic protease cascade. <i>Cell</i> 1997;91:479-89</p> <p>Hu Y, Benedict MA, Wu D, Inohara N, Nunez G. Bcl-XL interacts with Apaf-1 and inhibits Apaf-1-dependent caspase-9 activation. <i>Proc Natl Acad Sci USA</i> 1998;95:4386-91 (RD1299)</p> |

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