

ADAP (CT) antibody (pAb)

Rabbit Anti-Human/Mouse ADAP (CT)

Instruction Manual

| | |
|---------------------|--|
| Catalog Number | PK-AB718-4277 |
| Synonyms | ADAP Antibody: Adhesion and degranulation adapter protein, FYN binding protein, FYB-120, 130, SLAP-130 |
| Description | The adhesion and degranulation adaptor protein (ADAP) was initially identified as a molecular adapter that couples T cell receptor (TCR) stimulation to the avidity of integrins governing T cell adhesion. TCR stimulation promotes the formation of a multi-protein complex containing CARMA1, MALT1, and BCL-10, which through the association of ADAP, ultimately activates the NF-kappaB family of transcription factors. More recent experiments have shown that ADAP controls optimal T cell proliferation, cytokine production, and expression of the Bcl-2 family member Bcl-x(L), suggesting that ADAP regulates T cell activation by promoting antigen-dependent T cell-antigen presenting cell (APC) activation. At least three isoforms of ADAP are known to exist. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | ADAP antibody was raised in rabbits against a 20 amino acid peptide from near the carboxy terminus of human ADAP. |
| Purification Method | Affinity chromatography purified via peptide column. |
| Clone / IgG Subtype | Polyclonal antibody |
| Species Reactivity | Human, Mouse |
| Specificity | |
| Formulation | Antibody is supplied in PBS containing 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 | Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures. |
| Applications | E, WB, ICC, IF I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. ADAP antibody can be used for detection of ADAP by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL. |
| Images | Available upon request. |
| References | Griffiths EK, Krawczyk C, Kong YY, et al. Positive regulation of T cell activation and integrin adhesion by the adapter Fyb/Slap. <i>Science</i> 2001; 293:2260-3. Rawlings DJ, Sommer K, and Moreno-Garcia ME. The CARMA1 signalosome links the signalling machinery of adaptive and innate immunity in lymphocytes. <i>Nat. Rev. Immunol.</i> 2006; 6:799-812. Medeiros RB, Burbach BJ, Mueller KL, et al. Regulation of NF-kappaB activation in T cells via association of the adapter proteins ADAP and CARMA1. <i>Science</i> 2007; 316:754-8. Mueller KL, Thomas MS, Burbach BJ, et al. Adhesion and degranulation-promoting adapter protein (ADAP) positively regulates T cell sensitivity to antigen and T cell survival. <i>J. Immunol.</i> 2007; 179:3559-69. |
| Images | Available upon request. |
| Related Products | Cat.No. PK-AB718-4277P; ADAP Peptide Cat.No. PK-AB718-1204; K562 Cell Lysate |

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