

### Instruction Manual

|                                |  |
|--------------------------------|--|
| <b>Catalog Number</b>          | PK-AB718-2395  |
| <b>Synonyms</b>                | TACI Antibody:   |
| <b>Description</b>             | Members in the TNF superfamily regulate immune responses and induce apoptosis. Two novel members in the TNF family were recently identified and designated BAFF/BLyS/TALL-1/THANK/zTNF4 and April/TALL-2, respectively. BAFF was characterized as a B cell activator since it induced B cell proliferation and immunoglobulin secretion. April regulates immunological and non-immunological cell growth. Three receptors, BCMA, TACI, and BAFF-R, for BAFF and April were recently identified. TACI, like BCMA, binds BAFF and April. TACI and its ligands regulate humoral immune responses, activate NF-κB and c-jun N-terminal kinase, and are involved in B cell associated autoimmune diseases.                    |
| <b>Quantity</b>                | 100 µg   |
| <b>Source / Host</b>           | Rabbit   |
| <b>Subclass</b>                | Rabbit IgG   |
| <b>Immunogen</b>               | Rabbit polyclonal TACI antibody was raised against a synthetic peptide mapping at the amino terminus of human TACI.  |
| <b>Specific Species React.</b> | Human  |
| <b>Formulation</b>             | Antibody is supplied in PBS containing 0.02% sodium azide.   |
| <b>Stabilizer</b>              | sodium azide   |
| <b>Concentration</b>           | 1 mg/ml  |
| <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>Purification Method</b>     | Affinity chromatography purified via peptide column.   |
| <b>Applications</b>            | TACI polyclonal antibody can be used for detection of TACI by Western blot at 5 µg/mL. Antibody can also be used for immunocytochemistry starting at 2 µg/mL.  |
| <b>Storage</b>                 | E, WB, ICC   Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.<br>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.  |
| <b>References</b>              | Moore PA, Belvedere O, Orr A, et al. BLyS: member of the tumor necrosis factor family and B lymphocyte stimulator. <i>Science</i> 1999;285:260-3<br>Schneider P, MacKay F, Steiner V, et al. BAFF, a novel ligand of the tumor necrosis factor family, stimulates B cell growth. <i>J Exp Med</i> 1999;189:1747-56<br>Shu HB, Hu WH, Johnson H. TALL-1 is a novel member of the TNF family that is down-regulated by mitogens. <i>J Leukoc Biol</i> 1999;65:680-3<br>Mukhopadhyay A, Ni J, Zhai Y, Yu GL, Aggarwal BB. Identification and characterization of a novel cytokine, THANK, a TNF homologue that activates apoptosis, nuclear factor-κB, and c-Jun NH2-terminal kinase. <i>J Biol Chem</i> 1999 ;274:15978-81 |

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

# TACI antibody (pAb)

Rabbit Anti-Human TACI (Receptor for BAFF & APRIL)



| Related Products |   |
|------------------|---|
|                  | BAFF-R Antibody, Cat. No. PK-AB718-3097 |
|                  | TACI Peptide, Cat. No. PK-AB718-2395P   |
|                  | K562 Lysate, Cat. No. PK-AB718-1204     |
|                  | U937 Lysate, Cat. No. PK-AB718-1215     |

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

PromoCell GmbH  
Sickingenstr. 63/65  
69126 Heidelberg  
Germany

USA/Canada  
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)

02/2020