

### Instruction Manual

|                                |  |
|--------------------------------|--|
| <b>Catalog Number</b>          | PK-AB718-3023  |
| <b>Synonyms</b>                | Bcl-rambo Antibody; Bcl-2 like protein 13, Mil1  |
| <b>Description</b>             | Apoptosis plays a major role in normal organism development, tissue homeostasis, and removal of damaged cells. Disruption of this process has been implicated in a variety of diseases such as cancer. Members of the Bcl-2 family are known to be critical regulators of this process. These proteins are characterized by the presence of several conserved motifs termed Bcl-2 homology (BH) domains. A novel, widely expressed member termed Bcl-rambo was recently identified. This protein is localized to mitochondria in mammalian cells and its overexpression induces apoptosis which could be blocked by co-expression of inhibitor of apoptosis proteins (IAPs) such as XIAP, cIAP1, and cIAP2. Bcl-rambo shows overall homology to the anti-apoptotic members containing BH motifs, but unlike Bcl-2, the C-terminal membrane anchor of Bcl-rambo is preceded by a unique 250 amino acid insertion. This region by itself can induce apoptosis more efficiently than the Bcl-2 homology regions, suggesting that Bcl-rambo may be important other pro-apoptotic pathways. |
| <b>Quantity</b>                | 100 µg   |
| <b>Source / Host</b>           | Rabbit   |
| <b>Immunogen</b>               | Rabbit polyclonal Bcl-rambo antibody was raised against a 15 amino acid peptide from near the center of human Bcl-rambo (Genbank accession No. AAH07658).  |
| <b>Purification Method</b>     | Affinity chromatography purified via peptide column.   |
| <b>Clone / IgG Subtype</b>     | Polyclonal antibody  |
| <b>Species Reactivity</b>      | Human, Mouse, Rat  |
| <b>Specificity</b>             |  |
| <b>Formulation</b>             | Antibody is supplied in PBS containing 0.02% sodium azide.   |
| <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>Storage &amp; Stability</b> | 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>Applications</b>            | E, WB, IHC, IF <span style="float: right;">  Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.</span><br>Bcl-rambo antibody can be used for the detection of Bcl-rambo by Western blot at 2 and 4 µg/mL. Antibody can also be used for immunohistochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.   |
| <b>Images</b>                  | Available upon request.  |
| <b>References</b>              | Lockshin RA, Osborne B, and Zakeri Z. Cell death in the third millennium. <i>Cell Death Differ.</i> 2000; 7:2-7.<br>Cory S, Huang DCS, and Adams JM. The Bcl-2 family: roles in cell survival and oncogenesis. <i>Oncogene</i> 2003; 22:8590-607.<br>Heiser D, Labi V, Erlacher M, et al. The Bcl-2 protein family and its role in the development of neoplastic disease. <i>Exp. Gerontol.</i> 2004; 39:1125-35.<br>Kataoka T, Holler N, Michau O, et al. Bcl-rambo, a novel Bcl-2 homologue that induces apoptosis via its unique C-terminal extension. <i>J. Biol. Chem.</i> 2001; 276:19548-54.  |
| <b>Images</b>                  | Available upon request.  |
| <b>Related Products</b>        | Blocking Peptide, Cat. No. PK-AB718-3023P; K562 Cell Lysate, Cat. No. PK-AB718-1204<br>Bcl-2 Antibody (NT), Cat. No. PK-AB718-3335; cIAP Antibody (CT), Cat. No. PK-AB718-3325<br>XIAP Antibody (CT), Cat. No. PK-AB718-3331   |

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