



Anti- CARMA-3 / CARD-10, human (*Bimp1*)

CATALOG No.: 54984

BACKGROUND

CARMA proteins are scaffold molecules that belong to the membrane-associated guanylate kinase-like (MAGUK) and the caspase-associated recruitment domain (CARD) protein families. They play critical role in antigen receptor signaling pathways (1-4). CARMA1, CARMA2, and CARMA3 share homologous sequence and function but are found in different tissues. CARMA3 is most abundant in nonhematopoietic cells (5). CARMA3 protein physically associates with IKK complex and induces NF κ B activation via LPA signaling pathway (7). CARMA genes deactivation studies in mice resulted in reduced immunoglobulin levels and absence of TCR-induced NF κ B activation (4,6,7). This suggests that CARMA mutations may lead to adaptive and innate immunity malfunction.

SOURCE

Rabbit anti-human CARMA3 polyclonal antibody was raised against a synthetic peptide corresponding to the internal portion of human CARMA3. This sequence is also found in mouse, rat, chimpanzee, and monkey.

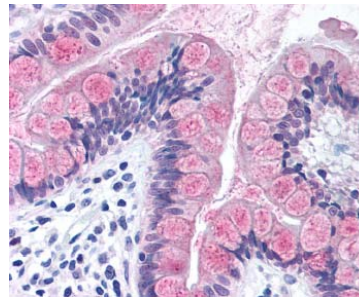
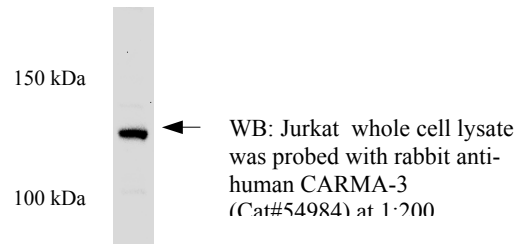
REACTIVITY

This epitope affinity purified rabbit polyclonal antibody reacts specifically with CARMA3. The antibody was evaluated by western blot, dot blot and ELISA. By western blot, an immunoreactive band around 130kDa was observed in NIH/3T3 and Jurkat whole cell lysates. In IHC, it showed moderate staining in a variety of formalin-fixed human tissues including heart, kidney, small and large intestine, testis, and skeletal muscle with minimal background staining. At least 22 different human formalin-fixed, paraffin archival tissues, and positive and negative tissues were scored and compared to the published literature on the expression and function of the gene.

APPLICATION

The antibody working concentrations for specific applications should be determined by the investigator. The following concentration ranges are recommended starting points for this product.

WB: 0.5 - 2 μ g/ml
IHC: 2.5-20 μ g/ml (10 μ g/ml-optimal)
IP*: 3.0-5.0 μ g/extract from 10^7 Cells
* Recommended but not tested.



Localization of the antibody as the precipitated red signal, with a hematoxylin purple nuclear counterstain in intestinal epithelium

This product is for in vitro research purposes only.

STORAGE

This polyclonal antibody is supplied as an epitope affinity purified rabbit IgG, 50 μ g in 250 μ l (0.2mg/ml) of 1X PBS (pH 7.4) containing 0.05% sodium azide. Store at 2-8 $^{\circ}$ C for up to one year from the date of shipment. Avoid repeated freezing thawing cycles.

REFERENCES

1. Gaide, O et al. (2002) Nature Immunology 3: 836-843
2. Hara, H. et al. (2004) JEM, 200(9): 1167-1177
3. Narayan, P. et al. (2006) Mol Cell Bio 26(6): 2327-2336
4. Egawa T. et al. (2003) Current Biology 13: 1252-1258
5. Grabiner, B. et al. (2007) Genes and Dev. 21: 984-996
6. Newton, K. et al. (2003) Current Biology 13: 1247-1251
7. Stilo R. et al. (2004) J. Biol. Chem. 279: 34323-34331