



Anti-ZAP-70

CATALOG NO.: 53958

BACKGROUND:

In humans, mutations in ZAP-70 gene results in a form of Severe Combined Immunodeficiency Syndrome (SCID). ZAP-70 expression also defines a subset of Chronic Lymphocytic Leukemia (CLL) in patients with unmutated Ig gene and poor clinical course.

SOURCE & REACTIVITY:

Mouse anti-ZAP-70 monoclonal antibody was raised against a recombinant protein encoding the amino acids 1-254 of human ZAP-70 with the immunoglobulin isotype being IgG2a. Anti-ZAP-70 reacts with ZAP-70 at the molecular weight of 70 kDa on western blot. Species reactivity includes human, while others are not tested.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

WB (Ab 1-2 µg/ml for 2hrs at room temperature (RT))

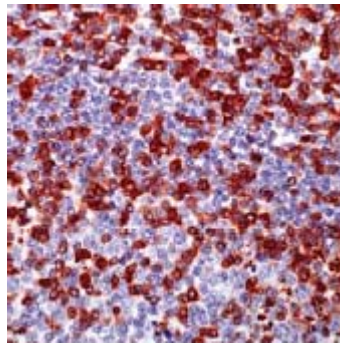
IHC (Formalin/paraffin)

Working dilution: 1:50 for 60 min at RT

Staining of formalin-fixed tissues requires boiling tissue sections in 10mM citrate buffer, pH 6.0 for 10 min followed by cooling at RT for 20 min.

Positive Control: T cells, NK cells, tonsil, jurkat cells.

Staining Pattern: Cytoplasmic (with occasional staining of macrophage lysozymes)



Human tonsil stained with anti-ZAP-70 (Cat. #53958)

This product is for in vitro research purposes only.

STORAGE:

0.25 ml antibody (at 200 µg/ml), purified from ascites fluid by Protein A, in 1X PBS (pH 7.4) with 0.2% BSA and 0.05% sodium azide. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles.