



Anti-Thrombospondin

CATALOG NO.: 53903

BACKGROUND:

Thrombospondin is a protein from platelet agranules. It is secreted at sites of platelet activation and aggregation and is involved in the processes of chemotaxis, adhesion, proliferation and differentiation of leukocytes, fibroblasts, smooth muscle and endothelial cells. This antibody inhibits TSP collagen interaction. Its binding to TSP is unaffected by glycosaminoglycans (e.g. hyaluronic acid, chondroitin sulfate, and heparin). It shows no cross reaction with fibronectin, fibrinogen, and von Willebrand factor. It shows a mild cross reactivity towards TSP 2.

SOURCE & REACTIVITY:

Mouse anti-thrombospondin monoclonal antibody was raised against a reduced and alkylated purified human TSP (fully denatured) from the supernatant of thrombin-activated platelets with the immunoglobulin isotype being IgG1 / κ . Anti-thrombospondin reacts with thrombospondin at the molecular weight of approximately 450 kDa as a non-reduced form and 170-180 kDa as a reduced form on western blot. Species reactivity includes human, mouse, rat, cow and pig, while others are not tested.

APPLICATION:

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

IP (Ab 2 μ g/mg protein lysate)

IFC

FC

WB (Ab 1-2 μ g/ml for 2hrs)

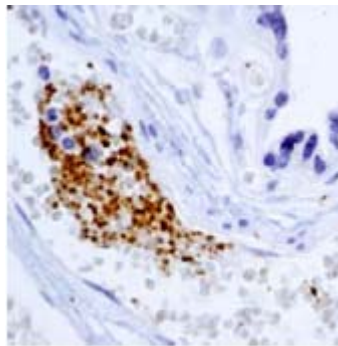
IHC (Formalin/paraffin)

Working dilution: 1:50 for 30 min at room temperature (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: Tonsil, placenta

Staining Pattern: Golgi complex, secretory granules, endoplasmic reticulum, extracellular matrix



Human placenta stained with anti-thrombospondin (Cat. #53903)

This product is for in vitro research purposes only.

STORAGE:

0.25 ml antibody (at 200 μ g/ml), purified from ascites fluid by Protein G, 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.