



Anti-PR3 (Proteinase3)

CATALOG NO.: 53791

BACKGROUND:

Proteinase 3 (PR3), a constituent of azurophil granules of neutrophils (polymorphonuclear cells, PMNs), is the target antigen for most anti-neutrophil cytoplasmic antibodies (c-ANCA) in Wegener's granulomatosis (WG).

SOURCE & REACTIVITY:

Anti-PR3 (proteinase3) mouse monoclonal antibody was raised against HMC-1/PR3 cell granule extract.

Molecular Weight: 32-40 kDa

Species reactivity includes human, while others remain untested.

APPLICATION:

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

IFC

Positive Control: HMC-1/PR3-S176 cells

This product is for in vitro research purposes only.

STORAGE:

This monoclonal antibody is supplied as 0.25 ml purified antibody from ascites fluid by protein G at 200 µg/ml in 10 mM PBS, pH 7.4, with 0.2% BSA and 0.05% sodium azide. Store at 2-8 °C for up to one year. Avoid repeated freezing and thawing.