

Purified Mouse Anti-human Myeloperoxidase Antibody *MPO421-8B2, monoclonal*

Catalog number: V1031805

Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Myeloperoxidase
Clone	MPO421-8B2

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED), ELISA

Applications

Myeloperoxidase (MPO) is a 84 kDa protein, located in the extracellular exosome, intracellular membrane-bounded organelle and nucleoplasm of cells. In Homo sapiens, myeloperoxidase suppresses apoptotic process, and moreover, recognizes heme, metal ion and chromatin. Myeloperoxidase is an integral part of organismal processes, in particular, removal of superoxide radicals, response to gold nanoparticle and response to lipopolysaccharide. It has been closely linked to critical functions like peroxidase activity. Mutations and abnormalities in myeloperoxidase have been closely linked to a number of diseases, in particular, myeloperoxidase deficiency (MPOD). Myeloperoxidase deficiency, an autosomal recessive inheritance disorder characterized by abnormality of blood and blood-forming tissues, abnormality of metabolism/homeostasis and abnormality of the immune system, has in specific been of interest to investigators.