

Purified Mouse Anti-Dextran Antibody
PolyclonalCatalog number: V102060
Unit size: 1.0 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	, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Mouse
Immunogen	Dextran

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	Flow Cytometry (FACS)

Applications

Dextran has shown use in a diverse array of applications, specifically in the field of biotechnological research and pharmacology. It has a variety of applications in the study of key cell-based receptors such as kinase receptors, especially in cell-lines like linking cells, other leukocytes and calcitriol-stimulated monocytes. Furthermore, dextran is vital to the examination of key biological proteins such as necroptotic proteins, VE-cadherin proteins and junctional protein. Upregulated activities, enzyme activities and prebiotic activity are, for instance, just a few of the activities that dextrans has helped scientists to better comprehend. It, in addition, has been widely used in the research of important cellular enzymes such as recombinant dextransucrase, co-immobilize dextransucrase and argininosuccinate synthetase. Dextran has a variety of applications in the research of essential biological pathways like gluconeogenetic pathway and Wnt pathways.