

**Purified Mouse Anti-human
immunodeficiency virus Capsid protein p24
(HIV-1) Antibody *ND1, monoclonal***

Catalog number: V103315
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 immunodeficiency virus
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Capsid protein p24 (HIV-1)
Clone	ND1

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, ICC, ELISA

Applications

Gag polyprotein is a protein with a molecular weight of 65 kDa, expressed in the cytoplasm, viral nucleocapsid and host multivesicular body of cells. In hiv-1s, gag polyprotein has been found to be involved in organismal processes, namely, viral budding via host ESCRT complex, RNA-dependent DNA biosynthetic process and nuclear transport. Gag polyprotein binds to RNA, zinc ion and N-myristoyl glycine. It has been associated with vital functions like structural molecule activity.