

**Purified Mouse Anti-human
immunodeficiency virus HIV protease
(HIV-1/HIV-2) Antibody *1696, monoclonal***

Catalog number: V103825
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	HIV protease (HIV-1/HIV-2)
Clone	1696

Biological Properties

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

Applications

Detects HIV-1/HIV-2 protease.