

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

Catalog number: V103820

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	Azide free Phosphate-buffered saline (PBS, pH 7.2) 0.2 µm filter sterilized, 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
Conjugate	Purified Azide Free

### Biological Properties

---

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

### Applications

---

Detects HIV-1/HIV-2 protease.