

**Purified Mouse Anti-human/mouse/pig
PSMA Antibody *GCP-04, monoclonal***

Catalog number: V1032055

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, mouse, pig, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PSMA
Clone	GCP-04

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, IHC(P), ICC

Applications

This is an antibody that targets PSMA.