

PE Mouse Anti-human/mouse/non-human primates PCNA Antibody *PC10, monoclonal*

Catalog number: V1031990

Unit size: 100 tests

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, non-human primates, drosophila, chicken, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PCNA
Clone	PC10
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

Applications

This is an antibody that targets PCNA.