

Purified Rabbit Anti-mycobacterium tuberculosis CFP-10 (M. tuberculosis) Antibody *PAb (490), monoclonal*

Catalog number: V103320

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	Mycobacterium tuberculosis
Class	Primary
Clonality	Monoclonal
Host	Rabbit
Immunogen	CFP-10 (M. tuberculosis)
Clone	PAb (490)

Biological Properties

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

Applications

ESAT-6-like protein is a protein with a molecular weight of 11 kDa.