

Purified Rabbit Anti-mycobacterium tuberculosis Tb7.7 (M. tuberculosis) Antibody *PAb (190), monoclonal*

Catalog number: V1032210
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	Tb7.7 (M. tuberculosis)
Clone	PAb (190)

Biological Properties

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

Applications

Antitoxin Rv2654c is a protein with a molecular weight of 9139 da. In m, antitoxin Rv2654c is a positive regulator of growth. (in M.smegmatis? based on what's on uniprot)