

Purified Rabbit Anti-mycobacterium tuberculosis ESAT-6 (M. tuberculosis) Antibody *PAb (491), monoclonal*

Catalog number: V103640
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	ESAT-6 (M. tuberculosis)
Clone	PAb (491)

Biological Properties

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

Applications

6 kDa early secretory antigenic target (alternatively called ESAT-6) is a 9904 da transmembrane protein that can be found in the cytoplasm, cell wall and integral component of membrane of cells. In m, 6 kDa early secretory antigenic target takes part in organismal processes, for instance, protein secretion by the type VII secretion system, suppression by symbiont of host T-cell mediated immune response and suppression by symbiont of host cell-mediated immune response. It binds to host cell surface. It is the subject of extensive research due to the fact that it is a component of the modulation by symbiont of host signal transduction pathway. It has been closely linked to critical functions like protein homodimerization activity.