

**Purified Rabbit Anti-mycobacterium tuberculosis GlcB (M. tuberculosis) Antibody \*PAb (192), monoclonal\***

Catalog number: V103740  
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	GlcB (M. tuberculosis)
Clone	PAb (192)

### Biological Properties

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

### Applications

Malate synthase G is a 80 kDa protein that can be found in the capsule, cytoplasm and plasma membrane of cells. In m, malate synthase G is thought to be essential to organismal processes, for instance, coenzyme A metabolic process, glyoxylate catabolic process and tricarboxylic acid cycle. Malate synthase G has been closely linked to vital functions such as malate synthase and protein homodimerization activity. It binds with host cell extracellular matrix, fibronectin and magnesium ion.