

**Purified Rabbit Anti-mycobacterium tuberculosis Acr1 (M. tuberculosis) Antibody \*PAb (185), monoclonal\***

Catalog number: V103015  
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	Acr1 (M. tuberculosis)
Clone	PAb (185)

### Biological Properties

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

### Applications

$\alpha$ -crystallin (Acr) is a 19 kDa protein that can be located in the extracellular region, cell wall and cytosol of cells. It is also known as Nox16 and HSP 16.3. Sequencing of HSP 16.3 has shown it contains a primary structural unit, the sHSP domain.  $\alpha$ -crystallin (Acr) aids in organismal processes, for example, cellular response to hypoxia and cellular response to starvation. It negatively regulates growth.