

**Purified Rabbit Anti-francisella tularensis
ssp. holarctica FTL0187 (F. tularensis ssp.
holarctica) Antibody *PAb (196),
monoclonal***

Catalog number: V103670
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 | Francisella tularensis ssp. holarctica |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Rabbit |
| Immunogen | FTL0187 (F. tularensis ssp. holarctica) |
| Clone | PAb (196) |

Biological Properties

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

Applications

Binds to Francisella tularensis subsp. holarctica cyclohexadienyl dehydratase.