

Purified Rabbit Anti-neisseria meningitidis Orf1 (N. meningitidis) Antibody *PAb (489), monoclonal*

Catalog number: V1031935

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 | Neisseria meningitidis |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Rabbit |
| Immunogen | Orf1 (N. meningitidis) |
| Clone | PAb (489) |

Biological Properties

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

Applications

Non-structural polyprotein pORF1 is a protein with a molecular weight of 216 kDa, found in the host cell cytoplasm of cells. In heV-1s, non-structural polyprotein pORF1 has been associated with essential functions like helicase, mRNA methyltransferase and RNA-directed 5'-3' RNA polymerase activity. Sequencing of non-structural polyprotein pORF1 has shown it contains 5 conserved structural units: alphavirus-like MT, RdRp catalytic, Macro, (+)RNA virus helicase C-terminal and (+)RNA virus helicase ATP-binding domain. Non-structural polyprotein pORF1 binds with identical protein, RNA and ATP. It has been found to be involved in organismal processes, namely, viral protein processing, DNA-templated transcription, and viral RNA genome replication.