

FITC Rabbit Anti-DYKDDDDK Antibody
Polyclonal

Catalog number: V102085

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 | , 0.2% (w/v) BSA |

Antibody Properties

| | |
|-----------|------------|
| Class | Primary |
| Clonality | Polyclonal |
| Host | Rabbit |
| Immunogen | DYKDDDDK |
| Conjugate | FITC |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions |
| Application | Flow Cytometry (FACS) |

Applications

Dykdddk peptide has exemplified use in a diverse array of applications, especially in the field of biomedicine and biochemistry. It is vital to the study of essential cell-based proteins like successfully detected FLAG-tagged proteins, especially in cell-lines like pluripotent stem cells.