

**Purified Rabbit Anti-DYKDDDDK Antibody**  
**\*Polyclonal\***Catalog number: V102075  
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   |

**Biological Properties**

---

|             |                                                                                                |
|-------------|------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with 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.