

**HRP Rabbit Anti-DYKDDDDK Antibody**  
**\*Polyclonal\***Catalog number: V102090  
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	HRP

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with HRP 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.