

Biotin Rabbit Anti-DYKDDDDK Antibody
PolyclonalCatalog number: V102080
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	Biotin

Biological Properties

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