

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
Formulation	HEPES – saline with 0.002% Thimerosal and stabilizer

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	DYKDDDDK

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with HRP under optimal conditions
Application	FACS

Spectral Properties

Conjugate	HRP
-----------	-----

Applications

Dykddddd 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.