

V5-tag Recombinant monoclonal antibody (SV5-P-K) (Human IgG1κ)

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS591-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Regulator of expression of viral proteins, Protein rev, GKPIP NPLLGLDST epitope tag, PK tag
Application	ELISA, IF, IHC (FS), IP, Radioimmunoassay, WB
Clone	SV5-P-K
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	Pk (V5) Epitope Tag (GKPIP NPLLGLDST).
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Virus
Specificity	Recognizes a 14 aa sequence (GKPIP NPLLGLDST) derived from the P and V proteins of the Paramyxovirus simian virus 5.
Technical Info / Product Notes	This is a chimeric human antibody, based on the original mouse IgG1 format, created for improved compatibility with existing reagents, assays and techniques.

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com