

PRODUCT INFORMATION



RSVF Protein Chimeric Monoclonal Antibody (Clone RSHZ19 (Felvizumab)) Item No. 37173

Overview and Properties

Contents:	This vial contains 200 µg of protein A-affinity purified monoclonal antibody
Synonyms:	Respiratory Syncytial Virus F Protein, Respiratory Syncytial Virus Fusion Protein, RSV Fusion Protein
Immunogen:	Recombinant human-targeted RSV F
Cross Reactivity:	(+) F protein
Species Reactivity:	(+) RSV; other species not tested
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 0.02% ProClin™ 300
Concentration:	1 mg/ml
Clone:	RSHZ19 (Felvizumab)
Host:	Chimeric Monoclonal Antibody
Isotype:	IgG1κ
Application:	ELISA; the optimal working concentration/dilution should be determined empirically.

Description

Respiratory syncytial virus (RSV) fusion (F) protein is a surface glycoprotein encoded by the *F* gene in RSV RNA.¹ It is synthesized as an inactive precursor protein, F_0 , that undergoes proteolytic cleavage to release the F_1 and F_2 subunits, which are joined together by two disulfide bonds.² Mature RSV F protein is composed of an N-terminal fusion peptide (FP), two heptad repeats (HRs), a transmembrane domain, and a cytoplasmic tail and assembles into homotrimers on the virus surface.¹ Upon insertion of the FP in the target cell membrane, the HRs form a six-helical bundle (6-HB) that enables RSV to fuse with the target cell. RSV F protein is highly conserved between RSV subtypes A and B with approximately 90% amino acid identities.³ RSV is the most common causative agent of pediatric lower respiratory tract infections.⁴ Cayman's RSV F Protein Chimeric Monoclonal Antibody (Clone RSHZ19 (Felvizumab)) was produced recombinantly from the original humanized RSHZ19 antibody sequence, substituting the mouse antigen-binding domain with the rabbit IgG1 antigen-binding domain, and can be used for ELISA and in functional assays. The RSHZ19 antibody was generated by fusing human IgG1 constant domains to the antigen-binding domain of a mouse anti-RSV F protein monoclonal antibody.⁵

References

1. Graham, B.S. and Anderson, L.J. Challenges and opportunities for respiratory syncytial virus vaccines. *Curr. Top. Microbiol. Immunol.* **372**, 391-404 (2013).
2. Day, N.D., Branigan, P.J., Liu, C., *et al.* Contribution of cysteine residues in the extracellular domain of the F protein of human respiratory syncytial virus to its function. *Viol. J.* **3**, 34 (2006).
3. Choi, S.-H., Park, K.S., and Kim, Y.-J. Analysis of respiratory syncytial virus fusion protein from clinical isolates of Korean children in palivizumab era, 2009-2015. *J. Infect. Chemother.* **25(7)**, 514-519 (2019).
4. Nair, H., Nokes, D.J., Gessner, B.D., *et al.* Global burden of acute lower respiratory infections due to respiratory syncytial virus in young children: A systematic review and meta-analysis. *Lancet* **375(9725)**, 1545-1455 (2010).
5. Tempest, P.R., Bremner, P., Lambert, M., *et al.* Reshaping a human monoclonal antibody to inhibit human respiratory syncytial virus infection in vivo. *Biotechnology (N.Y.)* **9(3)**, 266-271 (1991).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/07/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM