

CERTIFICATE of ANALYSIS



RSV F Protein Chimeric Monoclonal Antibody (Clone RSHZ19 (Felvizumab))

Item No. 37173 • Batch No. 0725740 • Expires 08NOV2025

Overview

Formulation:	200 µg of protein A-affinity purified monoclonal antibody PBS with 0.02% Proclin 300
Concentration:	1 mg/ml
Storage/Stability:	-20°C/1yr
Applications:	ELISA
Cross reactivity:	((+))F protein
Species reactivity:	((+))RSV

Reviewed and approved by: Emily Tucker

Approved date: 06NOV2024

A handwritten signature in black ink, appearing to read "Emily Tucker".

WARNING
THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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