

CERTIFICATE of ANALYSIS



SARS-CoV-2 Spike Glycoprotein Receptor Binding Domain L452R, T478K variant (rabbit IgG1 Fc-tagged)

Item No. 34837 • Batch No. 0638249 • Expires 30SEP2022

Overview

Formulation:	PBS, pH 7.4, with 5% mannitol, 5% D-(+)-trehalose, 0.01% Tween 20, and 10% glycerol
Concentration:	0.64 mg/mL
Storage/Stability:	-80°C/1yr
Purity:	≥90% estimated by SDS-PAGE
Association rate constant:	38,870.00
Dissociation rate constant (1/s):	0.00
Dissociation equilibrium constant:	7.93 nM
Maximum analyte response (RU):	245.60

Reviewed and approved by: Laura Kostrzewa

Approved date: 11FEB2022

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