

Certificate of Analysis

Product:	SeroFlash™ SARS-CoV-2 IgM/IgG Antibody Detection Kit
Catalog No.:	D-1001-025
Lot No.:	2004003
Description:	Rapid anti-SARS-CoV-2 IgG/IgM antibody test
Application:	For anti-SARS-CoV-2 IgG/IgM detection in human serum, plasma, whole blood, and fingertip blood
Test Method:	As described in the user guide
Size:	25 tests
Components:	Test cards, Dilution buffer
Storage:	Store at 2-30°C
Shelf Life:	Stable for 12 months from the date of shipment, when stored properly
Appearance:	Rectangular plastic white cassette measuring 70mm x 20mm x 5mm with centered rectangular window and off-centered round window
Cleanliness:	Reagents are clean and free from mold-releasing agents
Intended Use:	For Use by Healthcare Professionals or Research Scientists Only

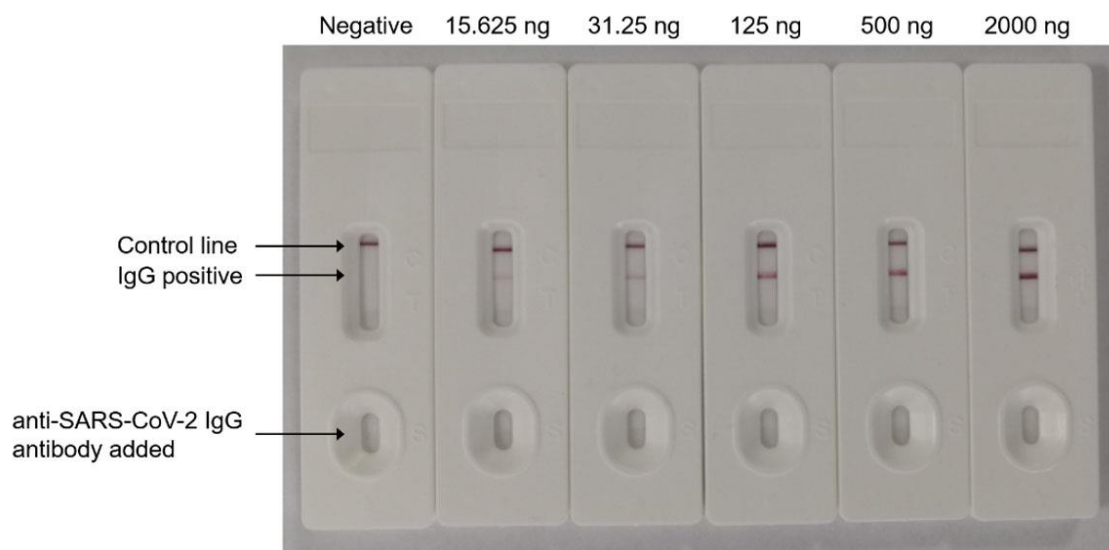


Figure 1. Detection of anti-SARS-CoV-2 IgG antibody at different concentrations with spiked humanized anti-SARS-CoV-2 IgG antibody.

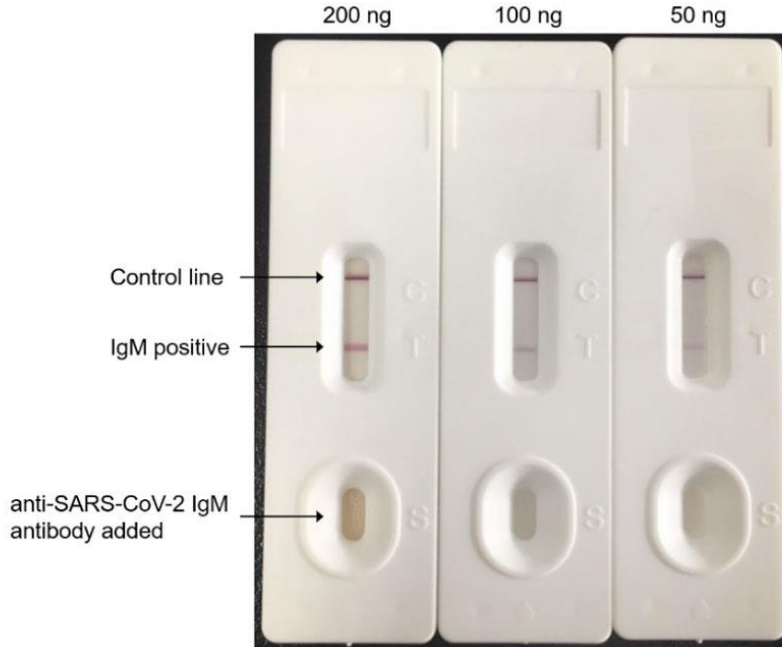


Figure 2. Detection of anti-SARS-CoV-2 IgM antibody at different concentrations with spiked humanized anti-SARS-CoV-2 IgM antibody.

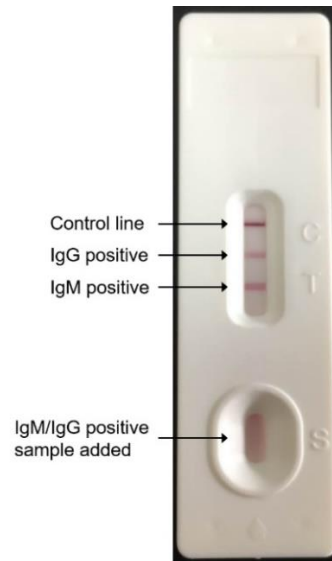


Figure 3. Detection of anti-SARS-CoV-2 IgG and IgM antibodies with clinically confirmed positive serum sample.

Approval: *Jin* Date: April 20, 2020

A signature on this CoA indicates that this product's lot meets quality specifications.