

ELISA PRODUCT INFORMATION & MANUAL

Human Respiratory Syncytial Virus Glycoprotein F - A2, (Prefusion) ELISA Kit (Colorimetric)

NBP3-26917

Sample Insert for reference use only

Enzyme-linked Immunosorbent Assay for quantitative detection. For research use only.

Not for diagnostic or therapeutic procedures.

PRINCIPLE OF THE ASSAY

The principle of this ELISA kit is based on the Sandwich Enzyme-Linked Immunosorbent Assay (ELISA). The 96-well strip plate is precoated with a monoclonal antibody specific for Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion). Standards and Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion) present in the sample is bound by the immobilized antibody. After washing, a horseradish peroxidase conjugated anti-Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion) monoclonal antibody is added, producing an antibody-antigen-antibody "sandwich complex". Following a wash to remove any unbound antibody-enzyme reagent, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion) bound in the initial step. The color development is stopped and the intensity of the color is measured at 450 nm.

KIT CONTENTS AND STORAGE

Please use the kit before the expiration date.

Components	Amt	Preparation instructions	storage
Pre-coated microplate	1plate (96 tests)	Take the microplate strips as needed, and put the unused strips back to the vacuum bag. It is best to vacuumize them.	Vacuum storage can store at 2-8°C until expiration date and opened package store at 2-8°C for one month.
Standard	1 bottle	Add 1 mL of 1 \times Dilution Buffer to the lyophilized standard bottle, briefly vortex to mix completely and prepare a standard stock solution.	The unopened lyophilized standard can be stored at 2-8°C until expiration date. Once the standard is redissolved, the liquid standard should be aliquoted and stored at -20°C to -80°C, please use it as soon as possible.
Detection Antibody	1 vial	Dilute at 1:660 with 1×dilution buffer for 10 minutes before use. Dilute fresh as needed.	
20× Dilution Buffer	1 bottle	If crystals have formed in the 20 \times concentrated solution, bring to room temperature and mix until dissolved. Dilute the 20 \times concentrated solution to	
20×Wash Buffer	1 bottle	$1 \times$ working solution with deionized water. For example, make 400 ml of $1 \times$ Wash Buffer by adding 20 mL of 20× Wash Buffer to 380 mL of deionized water. Dilute fresh as needed .	Primary liquid are stable at 2-8°C until expiration date. To be reconstituted, the working fluid is used within the working day and
Color Reagent A	1 bottle	Color Reagents A and B should be mixed together in equal volumes within 10 minutes before use. Take	discard. So dilute fresh as needed.
Color Reagent B	1 bottle	care not to contaminate the Color Reagent. If the mixed color reagent is blue, DO NOT USE .	
Stop Solution	1 bottle	Dilute acid. Use directly according to the use volume. Pay attention to safety when using.	

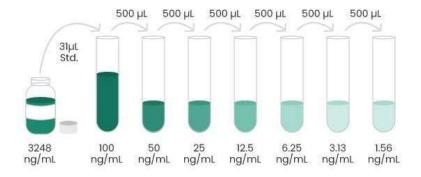
ASSAY PROCEDURE

1. Plate Set-up

- ➤ Bring all reagents to room temperature (22-28°C) equilibration (at least 30 minutes) before use. If crystals have formed in buffer solution, warm to room temperature and mix gently until the crystals have completely dissolved
- Determine the number of wells for the assay run. Remove unused microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and reseal.
- > Add 300μL1×Wash Buffer to each well and let stand for about 2 minutes. Aspirate or dump the liquid and pat dry on a paper towel, wash twice in this way.

2. Incubation with standard and samples [Volume: 100 μL Time: 1 hours]

Make standard curve: Prepare 1000 μL of the 100 pg/mL top standard by adding 31 μL of the standard stock solution in 969 μL of $1 \times \text{Dilution}$ Buffer. Perform six two-fold serial dilutions of the 100 pg/mL top standard in 6 separate tubes using 500 μL $1 \times \text{Dilution}$ Buffer as the diluents: after mixing the 100 pg/mL top standard, pipette 500 μL into the next tube, and so on. $1 \times \text{Dilution}$ Buffer serves as the zero standard (0 pg/mL). Ensures each assay has a standard curve. DO NOT USE the standard curve on other plates or other days.



- Add 100 μL standard and your test samples per well. Cover/seal the plate and incubate for 2 hours at room temperature.
- Add 300 µL 1×Wash Buffer to each well and let stand for about 2 minutes. Aspirate or dump the liquid and pat dry on a paper towel, wash wells 3 times in this way. Improper washes may lead to falsely elevated signals and poor reproducibility.

3. Incubation with Secondary Antibody [Volume: 100 μL Time: 0.5 hour]

- Add 100 μL of detection antibody working solution into each well, mix gently.
- Cover/seal the plate and incubate for 1 hour at room temperature.
- Removal the liquid in the wells and repeat the aspiration/wash as in Step 2.

4. Incubation with Substrate [Volume: 100 μL Time: about 20 minutes]

- Add 100 µL of Substrate Solution (the mixture of Color Reagents A and B) to each well, mix gently.
- Incubate for 20 minutes at room temperature. Protect from light. (According to the color of sample and the control antibody, the chromogenic time should be shortened or prolonged.)

5. Stop reaction

- Add 100 μL of Stop Solution to each well.
- > Tap gently the plate to ensure it is well mixed.

6. Absorbance Reading

> Read absorbance of the entire plate at 450nm wavelength within 10 minutes after adding the stop solution.

ASSAY PROCEDURE SUMMARY

1. Plate set-up Wash 2 times 2. Incubation with Add 100 µL standards or samples Incubate 2 hours, RT standard and samples Wash 3 times 3. Incubation with Add 100 µL Detection Antibody solution Secondary Antibody Incubate 1 hour, RT Wash 3 times 4. Incubation with Add 100 µL Substrate Solution Substrate Incubate 20 min, RT, in the dark 5. Stop reaction Add 100 µL Stop Solution 6. Absorbance Read absorbance at 450nm within 10 minutes Reading

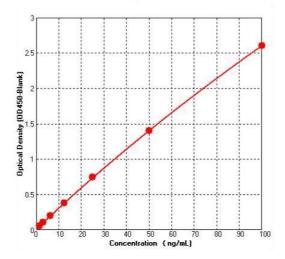
CALCULATION OF RESULTS

- 1. If samples generate values higher than the highest standard, dilute the samples and repeat the assay.
- 2. Calculate the mean absorbance for each standard and sample, subtract average zero standard optical density.
- 3. The data been calculated by 4-parameter logistics curve-fitting algorithm.

TYPICAL DATA

This standard curve is only for demonstration purposes. A standard curve should be run with each assay.

Concentration (ng/mL)	Zero standard subtracted OD-Blank
1.56	0.052
3.13	0.108
6.25	0.204
12.5	0.377
25	0.743
50	1.405
100	2.607



PERFORMANCE CHARACTERISTICS

Precision:

Intra-assay Precision (Precision within an assay) - Three samples of known concentration were tested twenty times on one plate to assess intra-assay precision.

Inter-assay Precision (Precision between assays) - Three samples of known concentration were tested in three separate assays to assess inter-assay precision.

	Intra -assay Precision			Inter -assay Precision		
Sample	1	2	3	1	2	3
N	20	20	20	3	3	3
Mean (pg/mL)	4	11	65	4	12	66
SD	0.24	0.76	1.97	0.05	0.30	2.62
CV (%)	6.6%	6.9%	3.1%	1.2%	2.6%	4.0%

Recovery: The recovery of Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion) spiked to different levels throughout the range of the assay in related matrices was evaluated.

Sample	Average % Recovery	Range
Cell culture supernates (n=3)	98	90 -108%

Linearity:

15	*	Cell culture supernates
1:2	recovery of detected	99%
1:4	recovery of detected	100%
1:8	recovery of detected	100%
1:16	recovery of detected	103%

Assay range: 1.56-100 pg/mL

Sensitivity: 0.1 pg/mL. Which was determined by adding two standard deviations to the mean optical density value of twenty zero standard replicates and calculating the corresponding concentration.

Specificity: This assay recognizes recombinant Human Respiratory Syncytial Virus Glycoprotein F - A2, (Pre-fusion).

PRECAUTIONS

- This kit is for research use only and is not for use in diagnostic or therapeutic procedures.
- 2. The kit should not be used beyond the expiration date.
- 3. Do not mix reagents from different lots.
- 4. The kit is designed and tested to detect the application which shown in the manual. The use of this kit for other purpose should be verified carefully by the end user.

SAFETY INSTRUCTIONS

- 1. The Stop Solution provided with this kit is an acid solution. Take care when using the reagent to avoid the risks
- 2. All biological materials should be handled and discarded as potentially hazardous following local laws and regulations.
- 3. Personal protective equipment such as lab coats, gloves, surgical masks and goggles are necessary in experiments for safety reasons.

TECHINICAL TIPS

- 1. Bring all reagents and samples to room temperature before use.
- 2. Samples should be thawed completely and mixed well prior to analysis. Avoid repeated freeze-thaw cycles of frozen samples.
- 3. Use a new disposable reagent reservoir and new disposable pipette tips for each transfer to avoid cross-contamination.
- 4. Read the absorbance of each well within 10 minutes after adding the stop solution.