



# SARS-COV2 CAPSID KITS

## PROTOCOL

### Part # 63ADK113PEG & 63ADK113PEH

**Test size#:** 500 tests (63ADK113PEG) and 10,000 tests (63ADK113PEH) - assay volume: 20  $\mu$ L

**Revision:** 02-Jan 2021

**Store at:** -60°C or below (63ADK113PEG); -60°C or below (63ADK113PEH)

**For research use only. Not for use in diagnostic procedures.**

### ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of SARS-CoV2 Capsid in cell lysates and offers a fast alternative to ELISA.

The detection principle of this kit is based on HTRF® technology (Homogeneous Time-Resolved Fluorescence). As shown in Figure 1, SARS-CoV2 Capsid is detected in a sandwich assay by using anti SARS-CoV2 capsid antibody labeled with Europium cryptate (donor), and anti SARS-CoV2 capsid antibody labeled with d2 (acceptor).

When the dyes are in close proximity, the excitation of the donor with a light source (laser or flash lamp) triggers a Fluorescence Resonance Energy Transfer (FRET) towards the acceptor, which in turn fluoresces at a specific wavelength (665 nm). Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the SARS-CoV2 Capsid concentration.

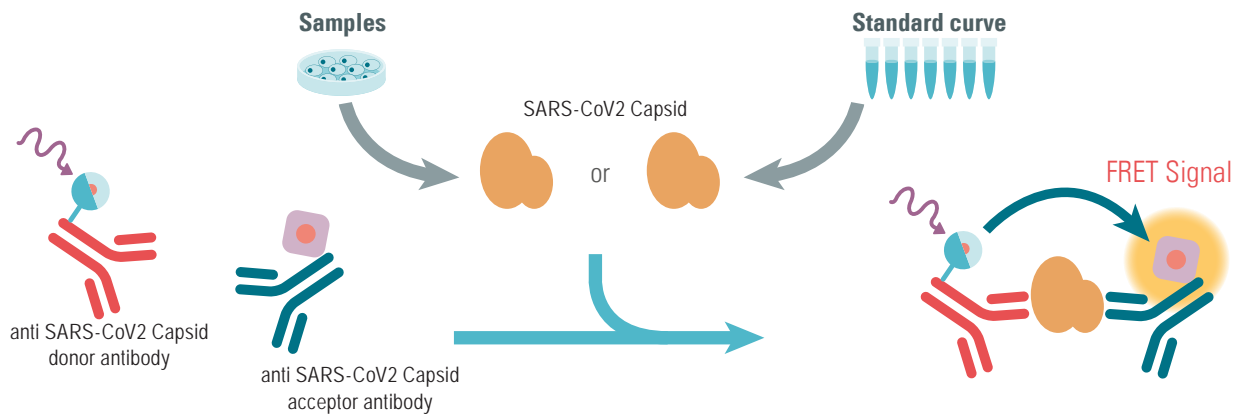
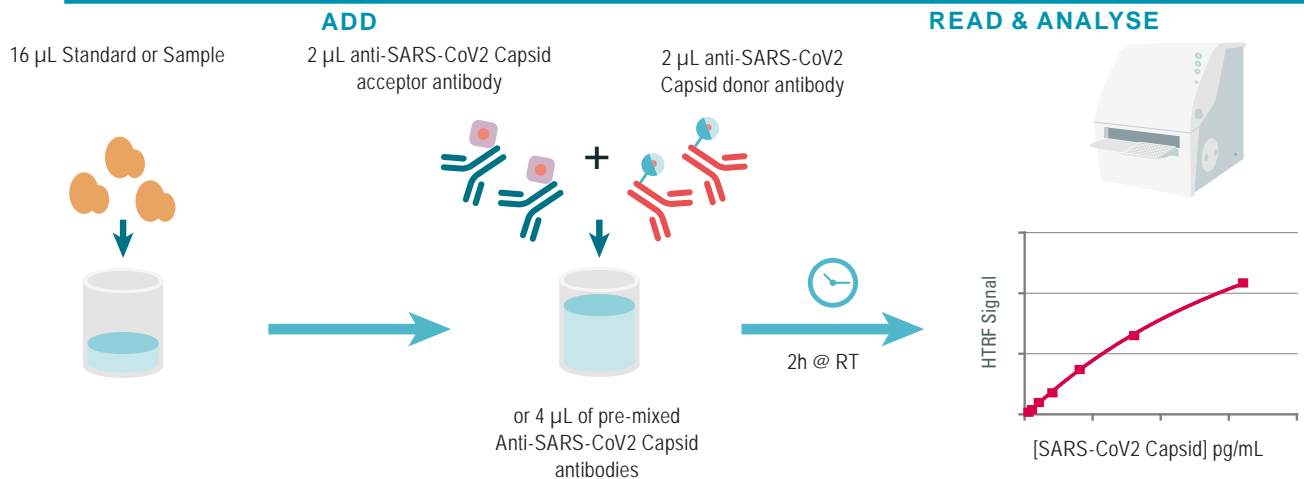


Figure 1: Principle of HTRF SARS-CoV2 Capsid sandwich assay.

### PROTOCOL AT A GLANCE



Make sure to use the set-up for Eu Cryptate.

**MATERIALS PROVIDED:**

<b>KIT COMPONENTS</b>	<b>500 TESTS * CAT # 63ADK113PEG</b>	<b>10,000 TESTS * CAT # 63ADK113PEH</b>
SARS-CoV2 Capsid Standard Frozen	1 vial - 50 µL 1.5 µg/mL	1 vial - 50 µL 1.5 µg/mL
SARS-CoV2 Capsid Eu Cryptate Antibody	1 vial - 20 µL Frozen - 50X	1 vial - 0.4 mL Frozen - 50X
SARS-CoV2 Capsid d2 Antibody	1 vial - 20 µL Frozen - 50X	1 vial - 0.4 mL Frozen - 50X
Lysis buffer #5 ** 4X	4 vials 2 mL	1 vial 130 mL
Detection buffer *** ready-to-use	1 vial 2 mL Detection Buffer	1 vial 50 mL Detection Buffer

\* When used as advised, the two available kit sizes will provide sufficient reagents for 500 tests and 10,000 tests respectively in 20 µL final volume..

Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

\*\* Medium like cell culture medium can be an alternative to the diluent.

\*\*\* The Detection buffer is used to prepare working solutions of acceptor and donor reagents.

**PURCHASE SEPARATELY:**

- HTRF®-Certified Reader. **Make sure the setup for Eu Cryptate is used.**

For a list of HTRF-compatible readers and set-up recommendations, please visit [www.cisbio.com/compatible-readers](http://www.cisbio.com/compatible-readers)

- Small volume (SV) detection microplates. For information about microplate recommendations, please visit our website at: [cisbio.com/microplates-recommendations](http://cisbio.com/microplates-recommendations)

**STORAGE AND STABILITY**

Store the kit at -60°C or below.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label. Detection buffer is shipped frozen, but can be stored at 2-8°C in your premises.



Reagents

If lyophilized, reconstituted reagents, antibodies, and standard stock solutions may be frozen and thawed only once. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below .

Volume of SARS CoV2 capsid standard aliquots should not be under 10 µL.

**REAGENT PREPARATION****BEFORE YOU BEGIN:**

- It is very important to prepare reagents in the specified buffers. The use of an incorrect lysis buffer may affect reagent stability and assay results.
- Thaw the frozen reagents at room temperature, allow them to warm up to room temperature for at least 30 mins before use
- Before use, allow Lysis buffer and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- SARS-CoV2 Capsid standards (for standard curve) must be prepared in lysis buffer or in the same medium as the samples.

**TAKE CARE TO PREPARE STOCK AND WORKING SOLUTIONS ACCORDING TO THE DIRECTIONS FOR THE KIT SIZE YOU HAVE PURCHASED.**

## TO PREPARE REAGENT STOCK SOLUTIONS:

500 TESTS KIT - 63ADK113PEG		10,000 TESTS KIT - 63ADK113PEH
Anti-SARS-CoV2 Capsid Eu Cryptate antibody		
Thaw the SARS-CoV2 Capsid Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.		Thaw the SARS-CoV2 Capsid Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.
Anti-SARS-CoV2 Capsid d2 antibody		
Thaw the SARS-CoV2 Capsid d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.		Thaw the SARS-CoV2 Capsid d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.
SARS-CoV2 Capsid Standard		
Thaw the SARS-CoV2 Capsid Standard solution (1.5 µg/mL) at RT. Mix gently. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solution into disposable plastic vials for storage at -60°C or below.		Thaw the SARS-CoV2 Capsid Standard solution (1.5 µg/mL) at RT. Mix gently. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solution into disposable plastic vials for storage at -60°C or below.
Lysis buffer		
Determine the amount of lysis buffer needed for the experiment. Each 96-well requires generally 50 µL of lysis buffer. Dilute 4-fold the 4 X lysis buffer with distilled water: homogenize the 4 X lysis buffer with a vortex and add 1 volume of stock solution in 3 volumes of distilled water (e.g., 1.25 mL of lysis buffer + 3.75 mL of distilled water). Mix gently after dilution. This 1 X lysis buffer is stable for 2 days at 2-8°C.		Determine the amount of lysis buffer needed for the experiment. Each 96-well requires generally 50 µL of lysis buffer. Dilute 4-fold the 4 X lysis buffer with distilled water: homogenize the 4 X lysis buffer with a vortex and add 1 volume of stock solution in 3 volumes of distilled water (e.g., 1.25 mL of lysis buffer + 3.75 mL of distilled water). Mix gently after dilution. This 1 X lysis buffer is stable for 2 days at 2-8°C.
Detection buffer		
The Detection buffer is ready-to-use.		The Detection buffer is ready-to-use.

## TO PREPARE ANTIBODY WORKING SOLUTIONS:

Each well requires 2 µL of SARS-CoV2 Capsid-Eu Cryptate Antibody and 2 µL of SARS-CoV2 Capsid-d2 Antibody.

Prepare the two antibody solutions in separate vials.

500 TESTS KIT - 63ADK113PEG		10,000 TESTS KIT - 63ADK113PEH
SARS-CoV2 Capsid Eu Cryptate antibody		
Dilute 50-fold the 50X stock solution (thawed reagent) of Anti-SARS-CoV2 capsid Eu Cryptate antibody with the Detection buffer: add 1 volume of Eu Cryptate antibody stock solution in 49 volumes of Detection buffer (e.g., 20 µL of reconstituted Eu Cryptate antibody stock solution + 980 µL of Detection Buffer).		Dilute 50-fold the 50X stock solution (thawed reagent) of Anti SARS-CoV2 capsid Eu Cryptate antibody with the Detection buffer: add 1 volume of Cryptate antibody stock solution in 49 volumes of Detection buffer (e.g., 0.4 mL of reconstituted Cryptate-antibody stock solution + 19.6 mL of Detection Buffer).
SARS-CoV2 Capsid d2 antibody		
Dilute 50-fold the 50X stock solution (thawed reagent) of Anti-SARS-CoV2 capsid d2 antibody with the Detection buffer: add 1 volume of d2 antibody stock solution in 49 volumes of Detection buffer (e.g., 20 µL of reconstituted d2 antibody stock solution + 980 µL of Detection Buffer).		Dilute 50-fold the 50X stock solution (thawed reagent) of Anti-SARS-CoV2 capsid d2 antibody with the Detection buffer: add 1 volume of d2 antibody stock solution in 49 volumes of Detection buffer (e.g., 0.4 mL of reconstituted d2 antibody stock solution + 19.6 mL of Detection Buffer).
Antibody mix		
It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).		It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).

## TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 16  $\mu\text{L}$  of standard.
- Dilute the standard stock solution serially with lysis buffer #5 (1X)
- In order to check for a potential interference effect from your own assay buffer when using the assay for the first time, we highly recommend the parallel preparation of a standard curve in your own supplemented cell culture medium and in lysis buffer #5 (1X).
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

A recommended standard dilution procedure is listed and illustrated below:

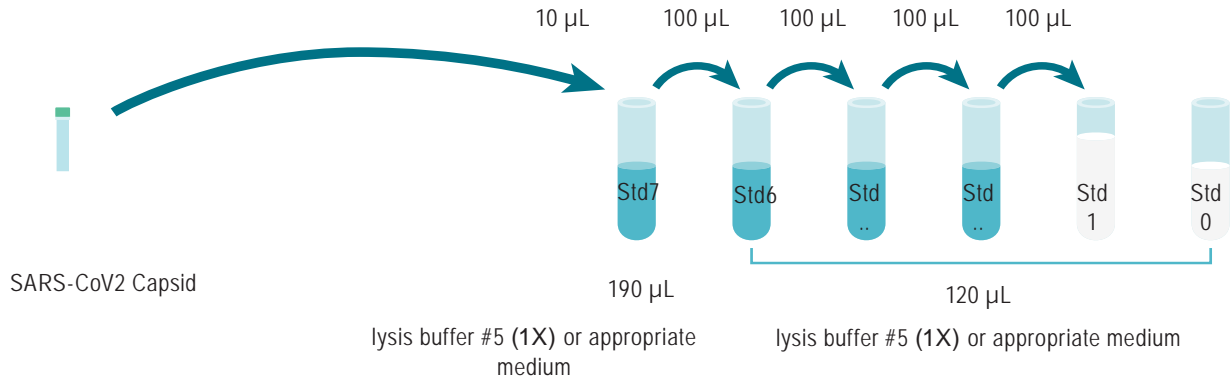
Dilute the standard stock solution 20-fold with lysis buffer; this yields the Standard Max solution (1.5  $\mu\text{g}/\text{mL}$ )

Dilute the standard stock solution 20-fold with lysis buffer #5 (1X) to prepare high standard (Std7): e.g. take 10  $\mu\text{L}$  of standard stock solution and add it to 190  $\mu\text{L}$  of lysis buffer #5 (1X). Mix gently.

Use the high standard (Std7) to prepare the standard curve using 1/2.2 serial dilutions as follows:

- Dispense 120  $\mu\text{L}$  of lysis buffer #5 (1X) in each vial from Std6 to Std 0.
- Add 100  $\mu\text{L}$  of standard to 120  $\mu\text{L}$  of lysis buffer #5 (1X), mix gently and repeat the 1/2.2 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 7 standards for the analyte. Std 0 (Negative control) is lysis buffer #5 (1X) or appropriate culture medium alone.








STANDARD	SERIAL DILUTIONS	SARS-COV2 CAPSID WORKING SOLUTIONS (PG/ML)
Standard Stock solution	Thawed stock solution	1 500 000
Standard 7	10 $\mu\text{L}$ Standard Stock Solution + 190 $\mu\text{L}$ lysis buffer 1X	75 000
Standard 6	100 $\mu\text{L}$ standard 7 + 120 $\mu\text{L}$ lysis buffer 1X	34 091
Standard 5	100 $\mu\text{L}$ standard 6 + 120 $\mu\text{L}$ lysis buffer 1X	15 496
Standard 4	100 $\mu\text{L}$ standard 5 + 120 $\mu\text{L}$ lysis buffer 1X	7 044
Standard 3	100 $\mu\text{L}$ standard 4 + 120 $\mu\text{L}$ lysis buffer 1X	3 202
Standard 2	100 $\mu\text{L}$ standard 3 + 120 $\mu\text{L}$ lysis buffer 1X	1 455
Standard 1	100 $\mu\text{L}$ standard 2 + 120 $\mu\text{L}$ lysis buffer 1X	661
Standard 0	120 $\mu\text{L}$ lysis buffer 1X	-

## TO PREPARE SAMPLES:

- Each well requires 16  $\mu$ L of sample.
- Just after their collection, put the samples at 4°C and test them immediately. For later use, samples should be dispensed into disposable plastic vials and stored at -60°C or below. Avoid multiple freeze/thaw cycles.
- Samples with a concentration above the highest standard (Std7) must be diluted lysis buffer #5 (1X)
- The assay can be run under a two-plate protocol, where cells are plated and stimulated in the same culture plate, then transferred to the assay plate for the HTRF® detection. This protocol enables the cells' viability and confluence to be monitored. It can also be further streamlined to a one-plate assay protocol where plating, stimulation and detection is performed in a single plate. For two-plate & one-plate assay protocols for suspension cells and adherent cells kept in medium for the lysis, we recommend to use the lysis buffer 4X (ready to use) For two-plate & one-plate assay protocols for adherent cells removing the medium for lysis, we recommend to use the lysis buffer 1X. Cell density, stimulation time, lysis step and other parameters related to the biology are cell-dependent and need to be optimized.
- To obtain additional information or support, please contact the HTRF technical support team at [cisbio.com/contact-us](https://cisbio.com/contact-us)

## ASSAY PROTOCOL

	Standard (Std 0 - Std7)	Samples
<b>Step 1</b> 	Dispense 16 $\mu$ L of each SARS-CoV2 Capsid standard (Std 0 - Std7) into each standard well	Dispense 16 $\mu$ L of each sample into each sample well
<b>Step 2</b> 	Add 2 $\mu$ L of SARS-CoV2 Capsid d2 antibody working solution to all wells	
<b>Step 3</b> 	Add 2 $\mu$ L of SARS-CoV2 Capsid Eu Cryptate antibody working solution to all wells	
<b>Step 4</b> 	Seal the plate and incubate 2h @ RT	
<b>Step 5</b> 	Remove the plate sealer and read on an HTRF® compatible reader	



## DATA REDUCTION

1. Calculate the ratio of the acceptor and donor emission signals for each individual well.

$$\text{Ratio} = \frac{\text{Signal 665 nm}}{\text{Signal 620 nm}} \times 10^4$$

2. Calculate the % CVs. The mean and standard deviation can then be worked out from ratio replicates.

$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

For more information about data reduction, please visit <http://www.cisbio.com/htrf-ratio-and-data-reduction>

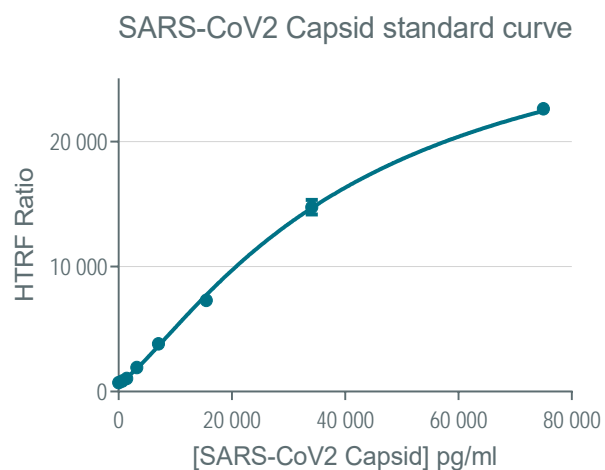
## RESULTS

This data must not be substituted for the data obtained in the laboratory and should be considered only as an example.

Results may vary from one HTRF® compatible reader to another.

Standard curve fitting with the 4 Parameter Logistic (4PL) model (with 1/Y<sup>2</sup> weighting):

	Ratio (1)	CV (2)
Standard 0 - Negative control	719	5%
Standard 1 - 661 pg/mL	835	4%
Standard 2 - 1,455 pg/mL	1050	9%
Standard 3 - 3,202 pg/mL	1917	8%
Standard 4 - 7,044 pg/mL	3817	0%
Standard 5 - 15,496 pg/mL	7308	3%
Standard 6 - 34,091 pg/mL	14759	4%
Standard 7 - 75,000 pg/mL	22637	2%



## ANALYTICAL CHARACTERISTICS

### ASSAY PERFORMANCES

Assay range (LOQ* to Std max)	307 - 75,000 pg/mL
Limit Of Detection (LOD)* = Mean Std 0 + 2 SD	165 pg/mL
Incubation time	2h at RT

\*The LOD and LOQ were calculated from data obtained in diluent with the PHERAstar FS reader (flash lamp excitation) after 2h incubation. These values may vary from one HTRF compatible reader to another.

**REACH European regulations and compliance** This product and/or some of its components include a Triton concentration of 0.1% or more and as such, it is concerned by the REACH European regulations. We recommend researchers using this product to act in compliance with REACH and in particular: to only use the product for in vitro research in appropriate and controlled premises by qualified researchers, ii) to ensure the collection and the treatment of subsequent waste, and iii) to make sure that the total amount of Triton handled does not exceed 1 ton per year.

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage. The use of the cell line will be done with appropriate safety and handling precautions to minimize health and environmental impact. Remaining disclaimer.

Copyright 2020 Cisbio Bioassays. All rights reserved. HTRF and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays.

#### FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880  
Japan +81-(0)43-306-8712 Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio