



# HUMAN AQ STING BINDING KITS

## PROTOCOL

**Part #** 64BDSTGQPEG & 64BDSTGQPEH

**Test size#:** 500 tests (64BDSTGQPEG), 10,000 tests (64BDSTGQPEH) - assay volume: 20  $\mu$ L

**Revision:** 02 / Oct. 2019

**Store at:** -60°C or below

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

### ASSAY PRINCIPLE

Cisbio Bioassays' Human AQ STING binding assay is only intended for quantitative measurement of human AQ STING ligand using HTRF<sup>®</sup> technology.

Human AQ STING ligand is detected in a competitive assay format using a specific Anti 6His antibody labeled with Terbium Cryptate (donor) which binds to human AQ STING protein 6His-tagged and AQ STING ligand labelled with d2 (acceptor). The detection principle is based on HTRF<sup>®</sup> technology. 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). Your compound competes with the AQ STING ligand labelled with d2, and thereby prevents FRET from occurring. The specific signal is inversely proportional to the compound concentration (Fig. 1).

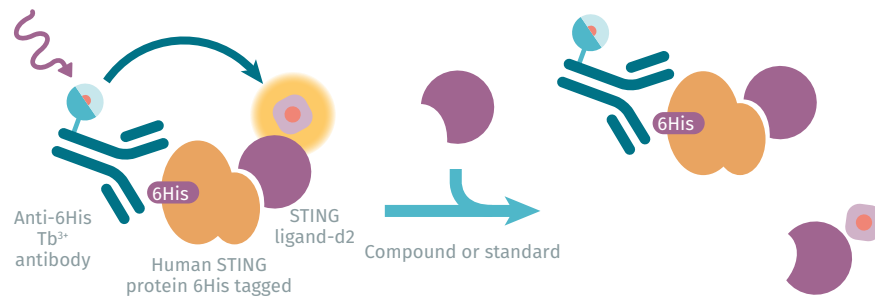
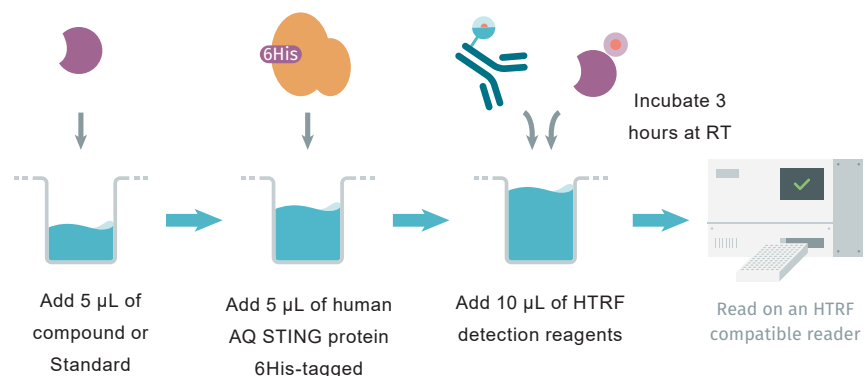


Figure 1: Principle of HTRF<sup>®</sup> Human AQ STING binding competitive assay.

### PROTOCOL AT A GLANCE



Make sure you use the appropriate setup for Tb<sup>3+</sup> Cryptate. For more information about setup and HTRF<sup>®</sup> compatible readers, please visit our website at: <http://www.cisbio.com/compatible-readers>

**MATERIALS PROVIDED:**

Kit components	500 tests * Cat # 64BDSTGQPEG	10,000 tests * Cat # 64BDSTGQPEH
Human AQ STING binding kit - Standard Frozen - 10X	1 vial - 50 $\mu$ L Ref 64BDSTGQCDA	2 vials - 50 $\mu$ L Ref 64BDSTGQCDA
Anti-6His-Tb <sup>3+</sup> Cryptate Antibody	1 vial - 50 $\mu$ L Frozen - 50 X	1 vial - 1 mL Frozen - 50 X
AQ STING ligand-d2	1 vial - 50 $\mu$ L Frozen - 50 X	1 vial - 1 mL Frozen - 50 X
human AQ STING protein 6His-tagged	2 vials - 25 $\mu$ L Frozen - 50 X	5 vials - 200 $\mu$ L Frozen - 50 X
Diluent #9 5X	3 vials 2 mL	1 vial 100 mL (62DL9DDC)
Detection buffer #12 5X	1 vial 2 mL	1 vial 50 mL

\* When used as advised, the two available kit sizes will provide sufficient reagents for 500 and 10,000 tests respectively in 20  $\mu$ L final. Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

**PURCHASE SEPARATELY:**

- HTRF<sup>®</sup> 96-well low volume plate Ref# 66PL96001 \*
- HTRF<sup>®</sup> 384-well low volume plate Ref 66PL384025 \*
- HTRF<sup>®</sup>-Certified Reader \*\*. Make sure the setup for Tb<sup>3+</sup> Cryptate is used.
- Use white plate only.

\* For HTRF microplate recommendations, please visit <http://www.cisbio.com/drug-discovery/htrf-microplate-recommendations>

\*\* For a list of HTRF-compatible readers and setup recommendations, please visit <http://www.cisbio.com/compatible-readers>

**STORAGE AND STABILITY**

Store the kit at -60°C or below. Under appropriate storage conditions, reagents are stable until the expiry date indicated on the label.



Reagents

Thaw and aliquot the protein on ice.

Once thawed, other solutions can be frozen 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 reagent aliquots should not be under 10  $\mu$ L.







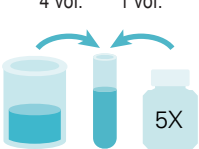
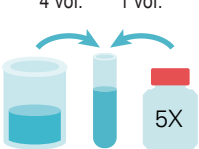
Thawed diluent and detection buffer can be stored at 2-8°C on your premises.

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

- It is very important to prepare reagents in the specified buffers. The use of an incorrect diluent may affect reagent stability and assay results.
- Thaw protein on ice, other reagents can be thawed at room temperature
- Before use, allow diluent and buffer to warm up at room temperature and homogenize them with a vortex.
- Human AQ STING binding kit - Standard (for standard curve) must be prepared in diluent.

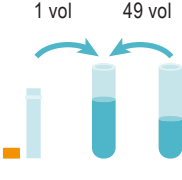
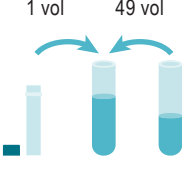


**TAKE CARE TO PREPARE STOCK AND WORKING SOLUTIONS ACCORDING TO THE DIRECTIONS FOR THE KIT SIZE YOU HAVE PURCHASED.**

## TO PREPARE REAGENT STOCK SOLUTIONS:

500 TESTS KIT - 64BDSTGQPEG			10,000 TESTS KIT - 64BDSTGQPEH
<b>Anti-6His- Cryptate antibody</b>			
Thaw the Anti-6His- Cryptate antibody. Centrifuge. This 50X stock solution can be frozen and stored at -60°C or below.			Thaw the Anti-6His- Cryptate antibody. Centrifuge. This 50X stock solution can be frozen and stored at -60°C or below.
<b>AQ STING ligand-d2</b>			
Thaw the AQ STING ligand-d2. Centrifuge. This 50X stock solution can be frozen and stored at -60°C or below.			Thaw the AQ STING ligand-d2. Centrifuge. This 50X stock solution can be frozen and stored at -60°C or below.
<b>Human AQ STING binding kit - Standard</b>			
Thaw the Human AQ STING binding kit - Standard. Centrifuge. This 10 X stock solution can be frozen and stored at -60°C or below.			Thaw the Human AQ STING binding kit - Standard. Centrifuge. This 10 X stock solution can be frozen and stored at -60°C or below.
<b>human AQ STING protein 6His-tagged</b>			
Thaw the human AQ STING protein 6His-tagged on ice. Centrifuge the vial. To avoid freeze/thaw cycles, it is recommended to aliquot the remainder of this 50X stock solution under 10 µL minimum in disposable plastic vials for storage at -60°C or below.			Thaw the human AQ STING protein 6His-tagged on ice. Centrifuge the vial. To avoid freeze/thaw cycles, it is recommended to aliquot the remainder of this 50X stock solution under 10 µL minimum in disposable plastic vials for storage at -60°C or below.
<b>Diluent</b>			
Dilute 5-fold the 5 X diluent #9 with distilled water: Homogenize the 5 X diluent #9 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water, e.g. 1 mL of diluent + 4 mL of distilled water. Mix gently after dilution.			Dilute 5-fold the 5 X diluent #9 with distilled water: Homogenize the 5 X diluent #9 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water, e.g. 10 mL of diluent + 40 mL of distilled water. Mix gently after dilution.
<b>Detection Buffer</b>			
Dilute 5-fold the 5 X detection buffer #12 with distilled water: Homogenize the 5 X detection buffer #12 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water, e.g. 1 mL of diluent + 4 mL of distilled water. Mix gently after dilution.			Dilute 5-fold the 5 X detection buffer #12 with distilled water: Homogenize the 5 X detection buffer #12 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water, e.g. 10 mL of diluent + 40 mL of distilled water. Mix gently after dilution.

**TO PREPARE WORKING SOLUTIONS:**

Each well requires 5  $\mu$ L of each reagent.  
Prepare in separate vials.

500 TESTS KIT - 64BDSTGQPEG	10,000 TESTS KIT - 64BDSTGQPEH	
<b>Anti-6His- Cryptate antibody</b>		
Dilute 50-fold the 50X stock solution (thawed reagent) of anti-6His cryptate antibody with detection buffer #12 (1X), eg 10 $\mu$ L of thawed cryptate-antibody stock solution + 490 $\mu$ L of detection buffer #12 (1X).		Dilute 50-fold the 50X stock solution (thawed reagent) of anti-6His cryptate antibody with detection buffer #12 (1X), eg 10 $\mu$ L of thawed cryptate-antibody stock solution + 490 $\mu$ L of detection buffer #12 (1X).
<b>AQ STING ligand-d2</b>		
Dilute 50-fold the 50X stock solution (thawed reagent) of AQ STING ligand-D2 with detection buffer #12 (1X), eg 10 $\mu$ L of thawed AQ STING ligand-D2 stock solution + 490 $\mu$ L of detection buffer #12 (1X).		Dilute 50-fold the 50X stock solution (thawed reagent) of AQ STING ligand-D2 with detection buffer #12 (1X), eg 10 $\mu$ L of thawed AQ STING ligand-D2 stock solution + 490 $\mu$ L of detection buffer #12 (1X).
<b>human AQ STING protein 6His-tagged</b>		
Dilute 50-fold the 50X stock solution (thawed reagent on ice) of human AQ STING 6His protein with detection buffer #12 (1X), eg 10 $\mu$ L of thawed protein stock solution + 490 $\mu$ L of detection buffer #12 (1X).		Dilute 50-fold the 50X stock solution (thawed reagent on ice) of human AQ STING 6His protein with detection buffer #12 (1X), eg 10 $\mu$ L of thawed protein stock solution + 490 $\mu$ L of detection buffer #12 (1X).
<b>HTRF reagents</b>		
It is possible to pre-mix the two ready-to-use solutions just prior to dispensing the reagents by adding 1 volume of AQ STING ligand-d2 solution to 1 volume of anti-6His cryptate antibody solution (e.g. 1 mL of AQ STING ligand-d2 + 1 mL of anti-6His cryptate antibody).		It is possible to pre-mix the two ready-to-use solutions just prior to dispensing the reagents by adding 1 volume of AQ STING ligand-d2 solution to 1 volume of anti-6His cryptate antibody solution (e.g. 1 mL of AQ STING ligand-d2 + 1 mL of anti-6His cryptate antibody).

### TO PREPARE WORKING STANDARD SOLUTIONS:

- Each well requires 5  $\mu\text{L}$  of standard.
- Dilute the standard stock solution serially with diluent #9 .
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

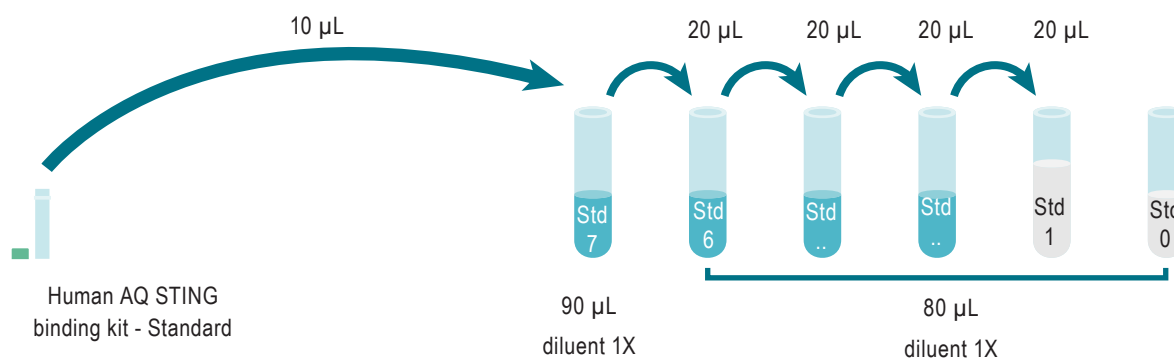
A recommended standard dilution procedure is listed and illustrated below:

Dilute the standard stock solution 10-fold with diluent to prepare high standard (Std 7): take 10  $\mu\text{L}$  of standard stock solution and add it to 90  $\mu\text{L}$  of diluent #9 (1X). Mix gently.

Use the high standard (Std 7) to prepare the standard curve using 1/5 serial dilutions as follows:

- Dispense 80  $\mu\text{L}$  of diluent into each vial from Std 6 to Std 0.
- Add 20  $\mu\text{L}$  of standard to 80  $\mu\text{L}$  of diluent, mix gently and repeat the 1/5 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 7 standards for the analyte. Std 0 (Positive control) is diluent alone.








STANDARD	SERIAL DILUTIONS	HUMAN AQ STING STANDARD WORKING SOLUTION (nM)	HUMAN AQ STING STANDARD FINAL CONCENTRATION (nM)
Standard Stock solution	Thawed stock solution	10 000	
Standard 7	10 $\mu\text{L}$ standard stock solution + 90 $\mu\text{L}$ Diluent	1 000	250
Standard 6	20 $\mu\text{L}$ standard 7 + 80 $\mu\text{L}$ Diluent	200	50
Standard 5	20 $\mu\text{L}$ standard 6 + 80 $\mu\text{L}$ Diluent	40	10
Standard 4	20 $\mu\text{L}$ standard 5 + 80 $\mu\text{L}$ Diluent	8	2
Standard 3	20 $\mu\text{L}$ standard 4 + 80 $\mu\text{L}$ Diluent	1.6	0.4
Standard 2	20 $\mu\text{L}$ standard 3 + 80 $\mu\text{L}$ Diluent	0.32	0.08
Standard 1	20 $\mu\text{L}$ standard 2 + 80 $\mu\text{L}$ Diluent	0.064	0.016
Standard 0	100 $\mu\text{L}$ Diluent	0	0

### TO PREPARE SAMPLES:

- Each well requires 5  $\mu\text{L}$  of compound.
- Dilute your compound in diluent #9 (1X).
- DMSO concentration must not exceed 2.5% final in the well (10% initial).

## ASSAY PROTOCOL

		<b>Negative control (or Cryptate control)</b>	<b>Standard (Std 0 - Std 7)</b>	<b>Compound</b>
<b>Step 1</b> 		Dispense 5 $\mu$ L of diluent into each negative control well.	Dispense 5 $\mu$ L of each Human AQ STING binding kit - Standard (Std 0 - Std 7) into each standard well.	Dispense 5 $\mu$ L of compound into each compound well.
<b>Step 2</b> 		Add 5 $\mu$ L of detection buffer to all wells	Add 5 $\mu$ L of human AQ STING protein 6His-tagged protein to all wells	
<b>Step 3</b> 	Add 10 $\mu$ L of AQ STING ligand-d2 and Anti 6His-Tb <sup>3+</sup> premixed working solution to all wells			
<b>Step 4</b> 	Seal the plate and incubate 3 hours at RT or at Over Night if necessary			
<b>Step 5</b> 	Remove the plate sealer and read on an HTRF <sup>®</sup> compatible reader			



## DATA REDUCTION & INTERPRETATION

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.

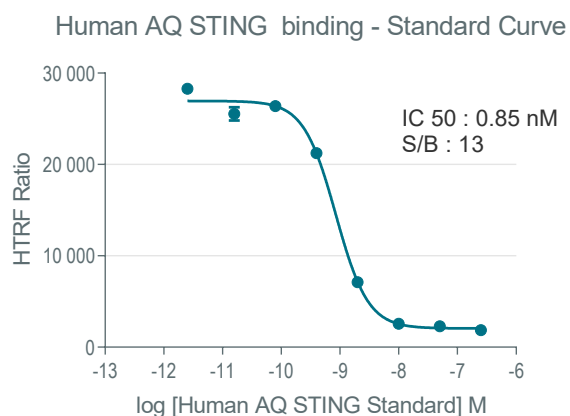
$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

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 (readouts on Pherastar FS with a flash lamp). Results may vary from one HTRF® compatible reader to another.

	Ratio <sup>(1)</sup>	CV <sup>(2)</sup>
Negative control	1844	1.5%
Std 0	28287	1.3%
Std 1 - 0.016 nM	25540	2.8%
Std 2 - 0.08 nM	26390	1.2%
Std 3 - 0.4 nM	21241	0.8%
Std 4 - 2 nM	7114	3.0%
Std 5 - 10 nM	2550	2.2%
Std 6 - 50 nM	2293	0.5%
Std 7 - 250 nM	1860	1.6%



## ANALYTICAL CHARACTERISTICS

Human AQ STING ligand-d2 Kd	14nM
Human AQ STING ligand-d2 concentration	14nM
Human AQ STING Standard IC50	0.85nM
Human AQ STING Standard Ki	0.42nM
Signal to background (S/B)	13

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.

Copyright 2018 Cisbio. All rights reserved.

### 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