

## Human IgG Lambda Kit 10,000 tests

[www.cisbio.com](http://www.cisbio.com)

### Product information

Document reference: 64LAMPEH-Rev02-Sept.2019

#### Packaging details:

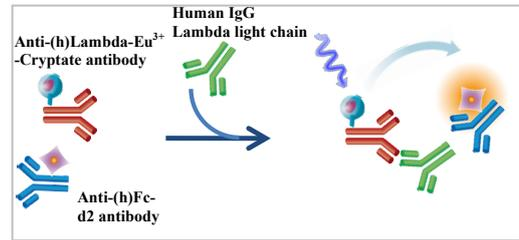
64LAMPEH	384-well low volume plate (20 µl) 10,000 tests
----------	---

### 1. ASSAY DESCRIPTION

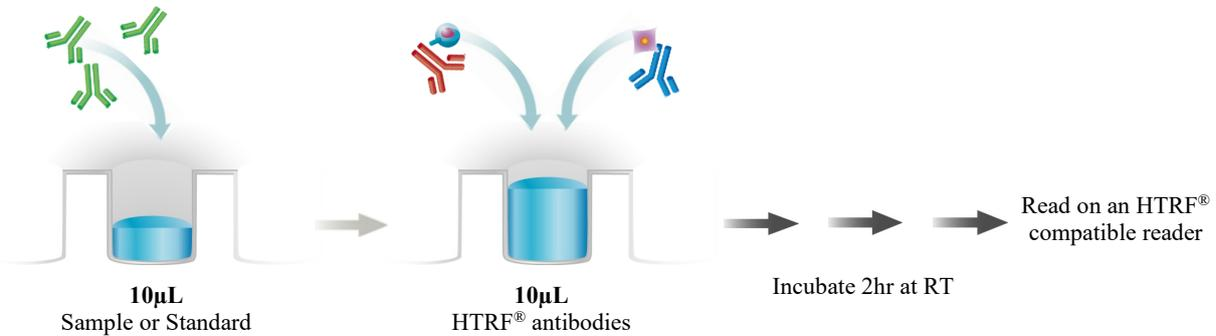
This assay is intended for the measurement of human (h) IgG Lambda light chain of all types of IgG (IgG1, IgG2, IgG3 and IgG4) using the HTRF<sup>®</sup> technology.

As shown here, (h)IgG Lambda light chain is detected in a sandwich assay format using 2 different specific antibodies. The anti-(h)-IgG Lambda antibody is labelled with Eu<sup>3+</sup>-Cryptate (donor) and the anti(h)-Fc antibody is labelled 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 (665nm). The two antibodies bind to the (h)IgG Lambda present in the sample, thereby generating FRET. The specific signal modulates positively in proportion to (h)IgG Lambda.



### 2. PROTOCOL AT A GLANCE



### 3. HTRF<sup>®</sup> REAGENTS

	Standard (h)IgGs	Anti-(h)Fc-d2 antibody	Anti-(h)IgG Lambda-Eu <sup>3+</sup> -Cryptate antibody	Diluent	Detection buffer #3
Stock solution	50 µL/vial 4 µg/mL	1000 µL/vial	1000 µL/vial	20 mL/vial	105 mL/vial
Storage	-20°C or below	-20°C or below	-20°C or below	4°C to -20°C*	4°C to -20°C*

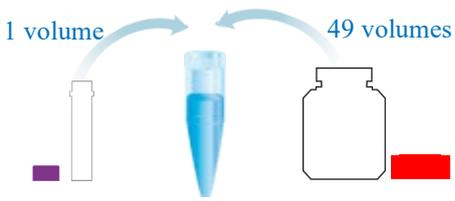
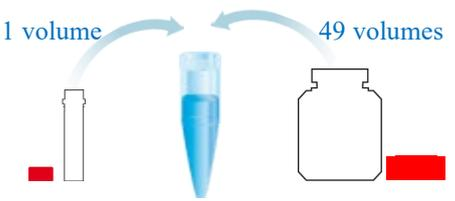
\* Diluent and Detection buffer are shipped frozen, but can then be stored at 2-8°C

## 4. REAGENT PREPARATION

Thaw all reagents at room temperature, allow them to warm up (caution: take buffers' thawing time into account).  
Prepare the working solutions from stock solutions (§3) by following the instructions below.

### 4.1. Preparation of antibody working solutions

Determine the amounts of antibodies needed for the experiment. Each well requires 5µL of each antibody.

Anti-(h)Fc-d2 antibody	Anti-(h)IgG Lambda-Eu <sup>3+</sup> -Cryptate antibody
	
Prepare a 50X diluted solution using the detection buffer #3: e.g. take 1000 µL of antibody stock solution and add it to 49 mL of detection buffer #3	Prepare a 50X diluted solution using the detection buffer #3: e.g. take 1000 µL of antibody stock solution and add it to 49 mL of detection buffer #3

Be careful, working solution preparation may differ between the 500 and the 10,000 data point kits.

### 4.2. Standard curve preparation

Determine how many standard levels and replicates to be tested. Each well requires 10µL of standard.

A whole IgG standard is provided with this kit.

For a more specific and quantitative calibration, we recommend the use of an appropriate IgG subtype: IgG1, IgG2, IgG3 or IgG4.

Standards	Working concentration (ng/mL)	Preparation
Std 9	200	20µL of Std stock solution + 380µL diluent
Std 8	100	100µl Std 9 + 100µl diluent
Std 7	50	100µl Std 8 + 100µl diluent
Std 6	25	100µl Std 7 + 100µl diluent
Std 5	12.5	100µl Std 6 + 100µl diluent
Std 4	6.25	100µl Std 5 + 100µl diluent
Std 3	3.1	100µl Std 4 + 100µl diluent
Std 2	1.6	100µl Std 3 + 100µl diluent
Std 1	0.8	100µl Std 2 + 100µl diluent
Std 0	0	100µl diluent

A recommended standard dilution procedure is listed below, and illustrated on the next page.

Dilute the standard stock solution 20-fold with diluent. This yields the high standard (Std 9) for the top of the curve (200ng/mL). In practice:

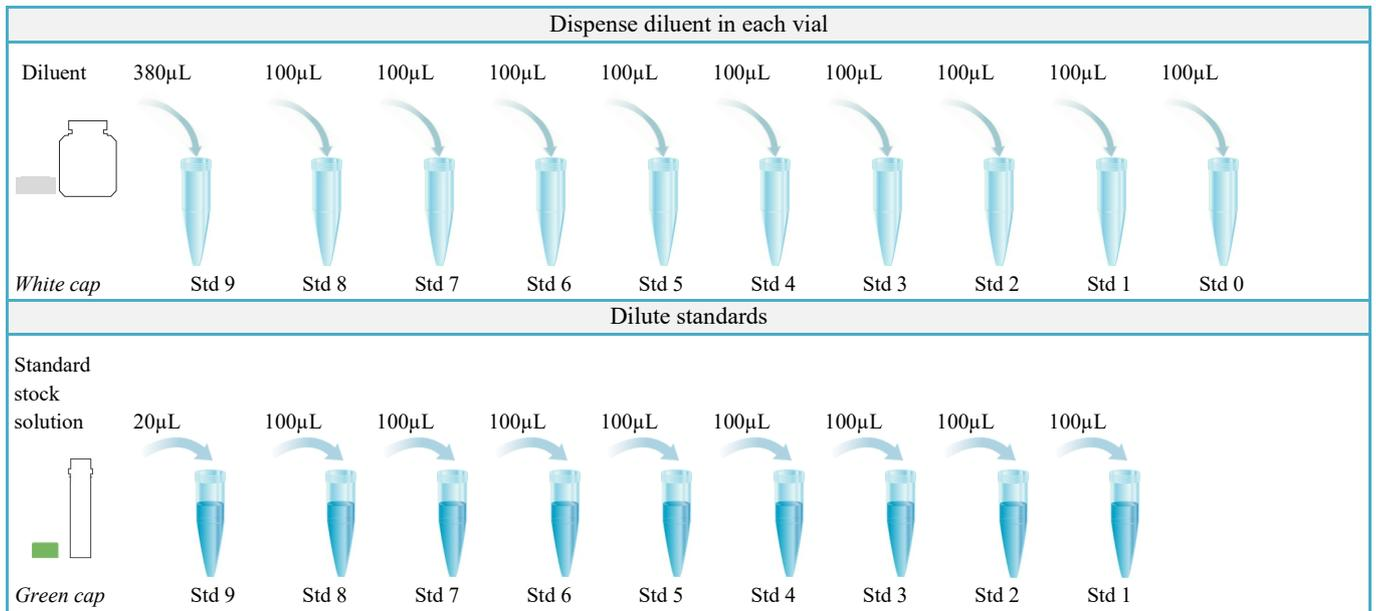
> e.g. take 20µL of the standard stock solution and add it to 380µL of diluent. Mix gently.

Use the high standard (Std 9) to prepare the standard curve using 1/2 serial dilutions as follows:

> Dispense 100µL of diluent in each vial from Std 8 to Std 1.

> Add 100µL of standard to 100µL of diluent, mix gently and repeat the 1/2 serial dilution to make standard solutions: 100, 50, 25, 12.5, 6.25, 3.1, 1.6, 0.8 ng/mL. This will create 9 standards for the analyte.

> Std 0 (negative control) is diluent alone.



#### Recommendations:

- HTRF<sup>®</sup> reagent concentrations have been set for optimal assay performances. Note that any dilution or improper use of the d2 and Cryptate antibodies will impair the assay's quality.
- For an accurate quantitative determination of sample, dilution must be carried out with the medium used for preparing the samples (i.e. diluent, culture medium or any other compatible medium).
- Standard and antibodies may be frozen and thawed once: to avoid freeze/thaw cycles it is recommended to dispense remaining stock solutions of standard and antibodies into disposable plastic vials for storage at -20°C or below.

To obtain additional information or support, please contact your technical support team ([htrfservices@cisbio.com](mailto:htrfservices@cisbio.com)).

## 5. ASSAY PROTOCOL

Dispense the reagents in the following order:



The 2 HTRF<sup>®</sup> antibodies can be pre-mix JUST PRIOR to dispensing: **DO NOT** store the pre-mix solution.

- ⇒ Cover the plate with a plate sealer
- ⇒ **Incubate at room temperature for 2 hours**
- ⇒ Remove the plate sealer and
- ⇒ Read the fluorescence emission at two different wavelengths (665nm and 620nm) on an HTRF<sup>®</sup> compatible reader (more information about compatible reader at [www.cisbio.com/readers](http://www.cisbio.com/readers))

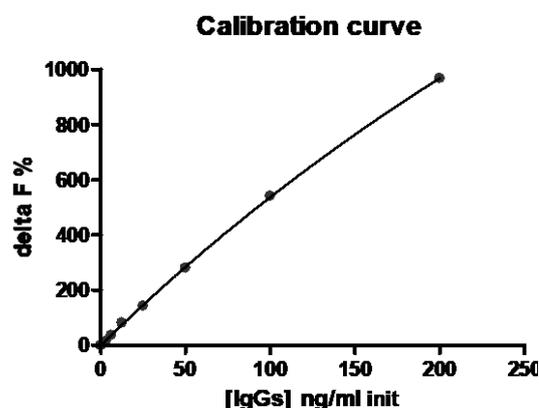
	Assay controls			Sample / Std
	Negative control	Cryptate control	Buffer control	
	<i>Used to calculate the delta F%</i>	<i>used to check the Cryptate signal at 620 nm</i>	<i>used to check background fluorescence</i>	
Sample / Std	-	-	-	10µL
Diluent	10µL	10µL	10µL	-
Anti-(h)Fc-d2 antibody	5µL	-	-	5µL
Anti-(h)IgG Lambda-Eu <sup>3+</sup> -Cryptate antibody	5µL	5µL		5µL
Detection buffer #3	-	5µL	10µL	-

## 6. DATA REDUCTION

This data must not be substituted for that obtained in the laboratory and should be considered only as an example (*readouts on PHERAstar<sup>plus</sup>*). Results may vary from one HTRF<sup>®</sup> compatible reader to another.

The assay standard curve is drawn up by plotting delta F% versus the analyte concentration:

Standard - ng/mL	Ratio (1)	CV % (2)	Delta F % (3)
Std 0 – Negative control	528	4.8%	0%
Std 1 – 0.8	532	3.6%	1%
Std 2 – 1.6	576	1.2%	9%
Std 3 – 3.1	603	2.1%	14%
Std 4 – 6.25	718	0.0%	36%
Std 5 – 12.5	956	2.4%	81%
Std 6 - 25	1286	2.2%	143%
Std 7 - 50	2007	2.9%	280%
Std 8 - 100	3390	0.3%	541%
Std 9 - 200	5645	1.6%	968%



Ratio (1)	$\frac{\text{Signal}_{665\text{nm}}}{\text{Signal}_{620\text{nm}}} \times 10^4$	Ratio must be calculated for each individual well.
CV % (2)	$\frac{\text{Standard deviation}}{\text{Mean ratio}} \times 100$	The mean and standard deviation can then be worked out from ratio replicates.
Delta F % (3)	$\frac{\text{Ratio}_{\text{standard or sample}} - \text{Ratio}_{\text{Negative control}}}{\text{Ratio}_{\text{Negative control}}} \times 100$	Reflects the signal to background of the assay. The negative control plays the role of an internal assay control.

For more information about data reduction, please visit our website at: [www.cisbio.com/data-reduction](http://www.cisbio.com/data-reduction)

## 7. ASSAY CHARACTERISTICS

### 7.1. Cross-reactivity

	Cross-reactivity %
Human Lambda	100
Human Kappa	0
Mouse Lambda	0
Human IgM	0

### 7.2. Detection limit

Human Lambda (IgG1) = 0.4ng/mL

To obtain additional information or support, please contact your technical support team ([htrfservices@cisbio.com](mailto:htrfservices@cisbio.com)).