

## Human IgG Lambda Kit 500 tests

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

### Product information

Document reference: 64LAMPEG-Rev02-Sept.2019

#### Packaging details:

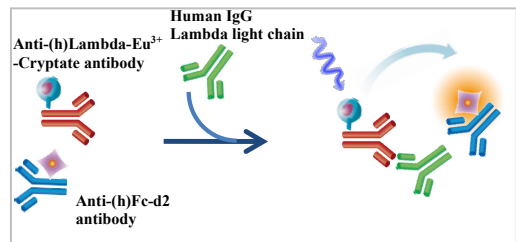
|          |  |
|----------|--|
| 64LAMPEG | 384-well low volume plate (20 µl)<br>500 tests |
|----------|--|

### 1. ASSAY DESCRIPTION

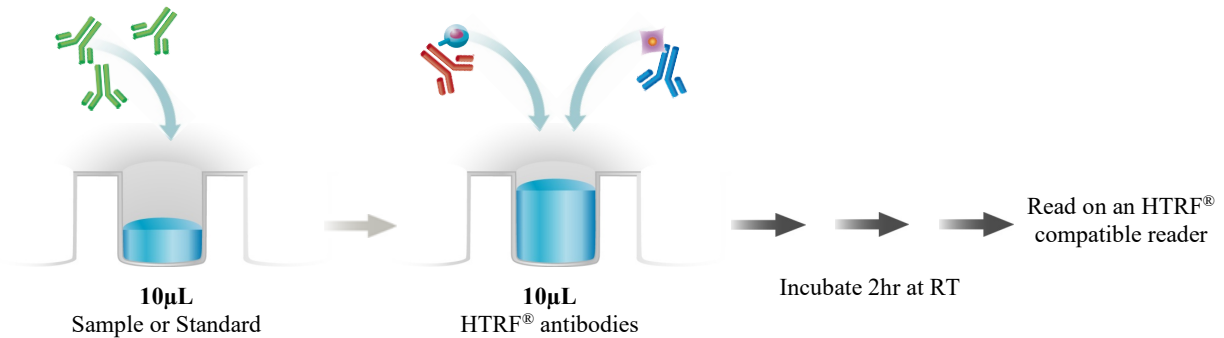
This assay is intended for the measurement of human (h) IgG Lambda light chain of all types of IgG (IgG1, IgG2, IgG3 and IgG4) using the HTRF® technology.

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® 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<br>4 µg/mL | 50 µL/vial             | 50 µL/vial   | 20 mL/vial    | 7 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*       |

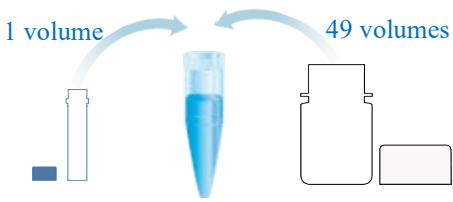
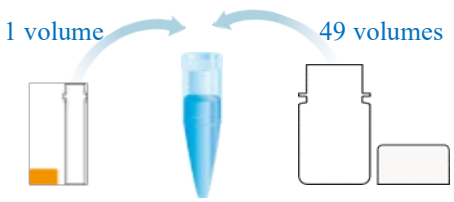
\* Diluent and Detection buffer are shipped frozen, but can then be stored at 2-8°C

## 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:<br>e.g. take 50 µL of antibody stock solution and add it to 2450 µL of detection buffer #3 | Prepare a 50X diluted solution using the detection buffer #3:<br>e.g. take 50 µL of antibody stock solution and add it to 2450 µL 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:

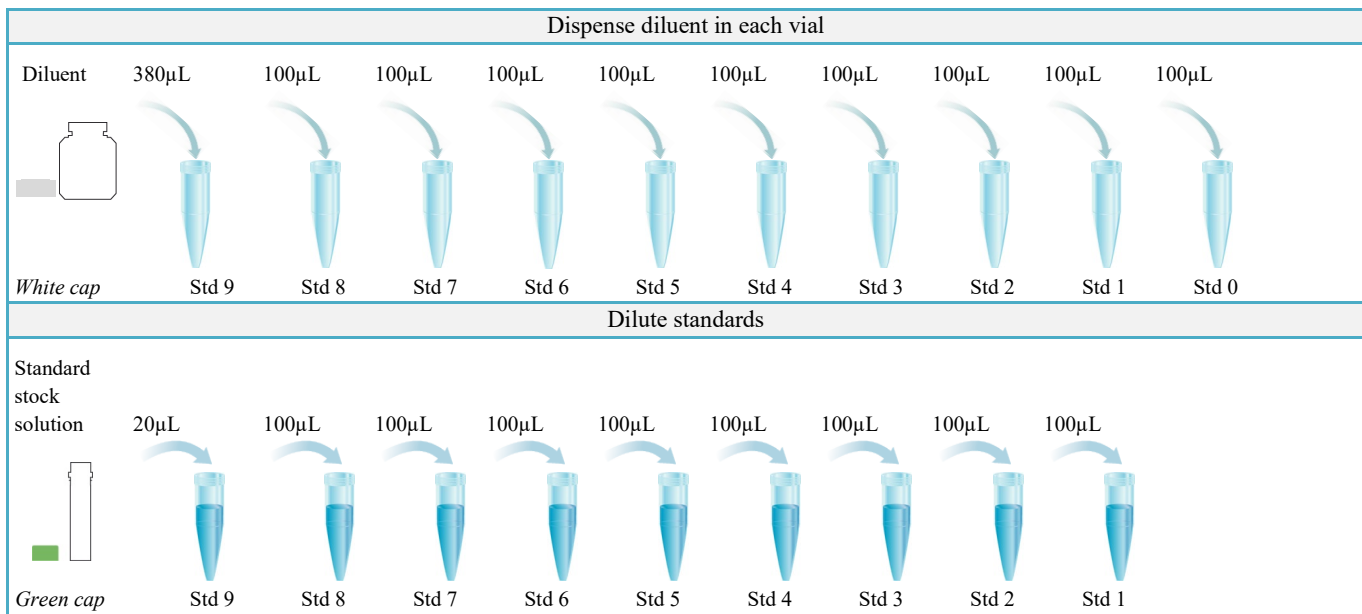
> e.g. take 20µL of the standard stock solution and add it to 380µL of diluent. Mix gently.

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.



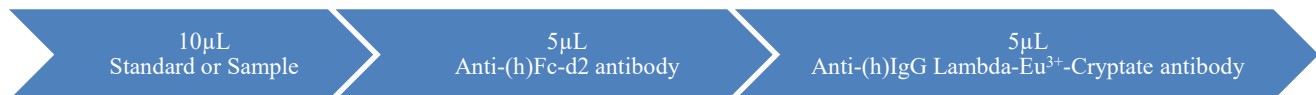
#### Recommendations:

- 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))

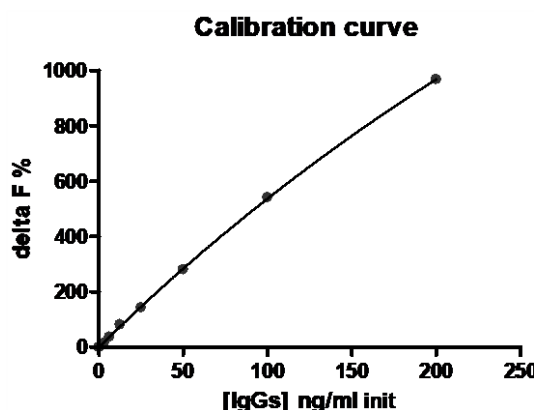
|  | Assay controls                        |  |  | Sample / Std |
|--|---------------------------------------|--|--|--------------|
|  | Negative control                      | Cryptate control                                   | Buffer control                               |              |
|  | <i>Used 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)).