

# MANUAL

Technology: HTRF®

## HTRF Human IgG Kappa kit

Part number	64KAPPEG	64KAPPEH
Test size	500 tests	10,000 tests

Storage:  $\leq -20^{\circ}\text{C}$

Version: 04

Date: January 2024

## ASSAY PRINCIPLE

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

As shown here, (h)IgG Kappa light chain is detected in a sandwich assay format using 2 different specific antibodies. The anti-(h)-IgG Kappa antibody is labelled with d2 (acceptor) and the anti(h)-Fc antibody is labelled with Eu3+-Cryptate (donor).

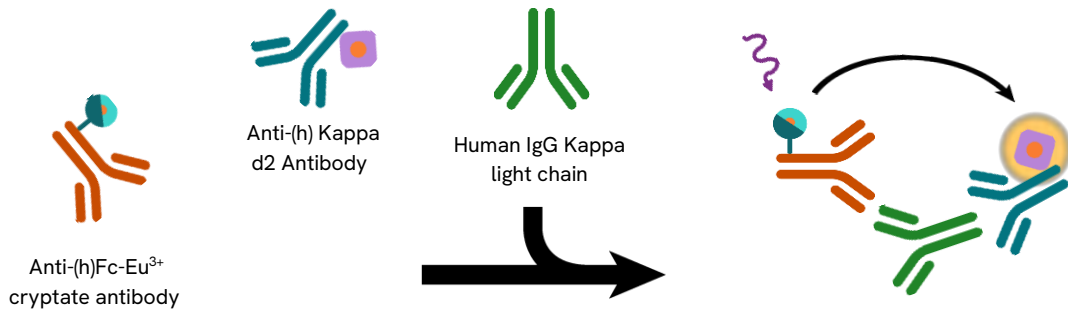
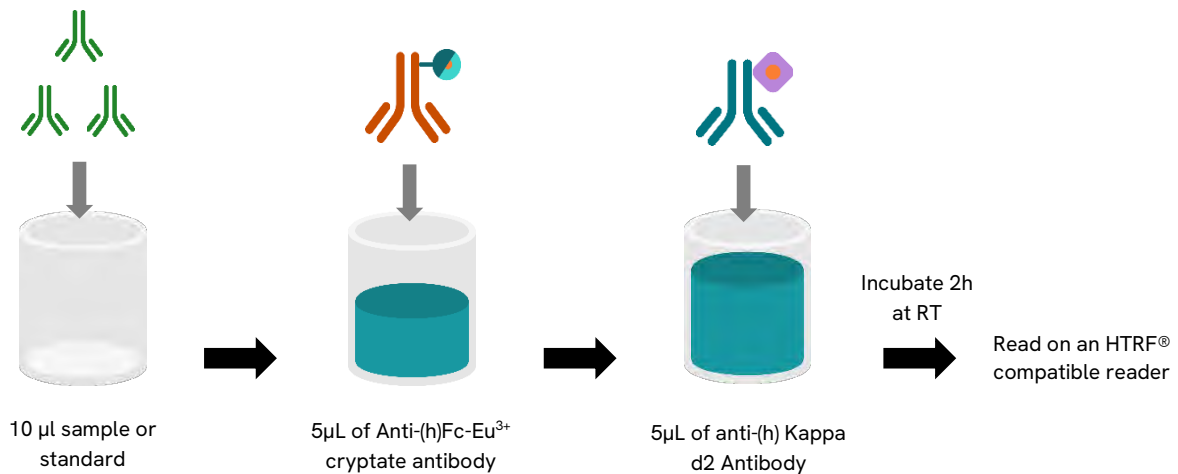












Figure 1: Principle of the assay.

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 Kappa present in the sample, thereby generating FRET. The specific signal modulates positively in proportion to (h)IgG Kappa.

## PROTOCOL AT A GLANCE



## MATERIAL PROVIDED

KIT COMPONENTS	STORAGE	500 TESTS			10,000 TESTS		
Standard (h)IgGs	≤-20°C		green cap	50 µl/vial 4 µg/mL		green cap	50 µl/vial 4 µg/mL
Anti-(h) IgG Kappa d2 antibody	≤-20°C		blue cap	50 µl/vial		purple cap	1,000 µl/vial
Anti-(h) Fc Eu <sup>3+</sup> Cryptate antibody	≤-20°C		orange cap	50 µl/vial		orange cap	1,000 µl/vial
Diluent	4°C to -20°C*		transparent cap	20 ml/vial		transparent cap	20 ml/vial
Detection buffer #3	4°C to -20°C*		red cap	7 ml/vial		red cap	105 ml/vial

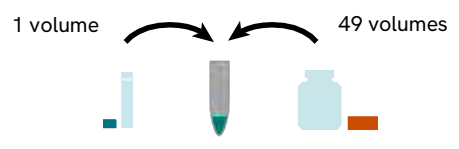
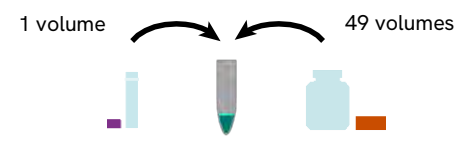
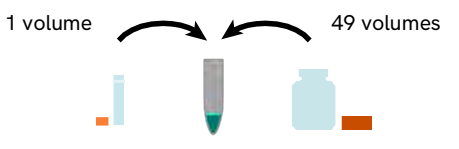
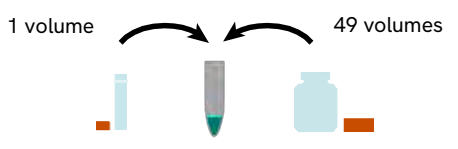
\* Diluent and Detection buffer are shipped frozen, but can then be stored at 2-8°C

## 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 by following the instructions below.

### Preparation of antibody working solutions

Determine the amounts of each detection reagent needed for the experiment. Each well requires 5µL of each detection reagent.

500 TESTS KIT	10,000 TESTS KIT
<b>Anti-(h) IgG Kappa-d2 antibody</b>	
	
Prepare a 50X diluted solution using the detection buffer#3: e.g. take 50 µL of detection reagent stock solution and add it to 2450 µL of detection buffer#3.	Prepare a 50X diluted solution using the detection buffer#3: e.g. take 1 mL of detection reagent stock solution and add it to 49 mL of detection buffer#3.
<b>Anti-(h) Fc-Eu<sup>3+</sup>-Cryptate antibody</b>	
	
Prepare a 50X diluted solution using the detection buffer#3: e.g. take 50 µL of detection reagent stock solution and add it to 2450 µL of detection buffer#3.	Prepare a 50X diluted solution using the detection buffer#3: e.g. take 1 mL of detection reagent stock solution and add it to 49 mL of detection buffer#3.

## Standard curve preparation

Determine how many standard levels and replicates to be tested. Each well requires 10  $\mu\text{L}$  of standard.

A whole IgG standard is provided with this kit.

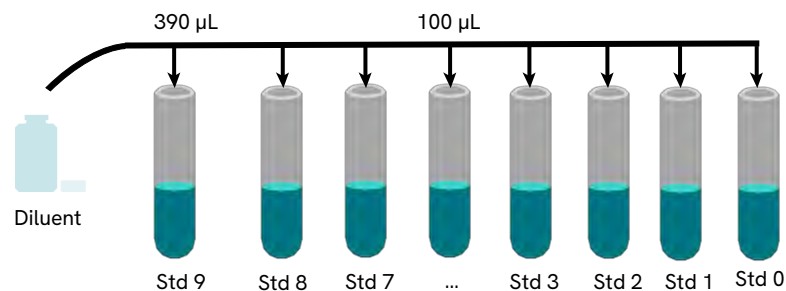
For a more specific and quantitative calibration, we recommend the use of an appropriate IgG subtype: IgG1, IgG2, IgG3 or IgG4.

STANDARD	PREPARATION	WORKING CONCENTRATION (ng/mL)
Standard 9	10 $\mu\text{L}$ of standard stock solution + 390 $\mu\text{L}$ diluent	100
Standard 8	100 $\mu\text{L}$ Std 9 + 100 $\mu\text{L}$ diluent	50
Standard 7	100 $\mu\text{L}$ Std 8 + 100 $\mu\text{L}$ diluent	25
Standard 6	100 $\mu\text{L}$ Std 7 + 100 $\mu\text{L}$ diluent	12.5
Standard 5	100 $\mu\text{L}$ Std 6 + 100 $\mu\text{L}$ diluent	6.25
Standard 4	100 $\mu\text{L}$ Std 5 + 100 $\mu\text{L}$ diluent	3.1
Standard 3	100 $\mu\text{L}$ Std 4 + 100 $\mu\text{L}$ diluent	1.6
Standard 2	100 $\mu\text{L}$ Std 3 + 100 $\mu\text{L}$ diluent	0.8
Standard 1	100 $\mu\text{L}$ Std 2 + 100 $\mu\text{L}$ diluent	0.4
Standard 0	100 $\mu\text{L}$ diluent	0

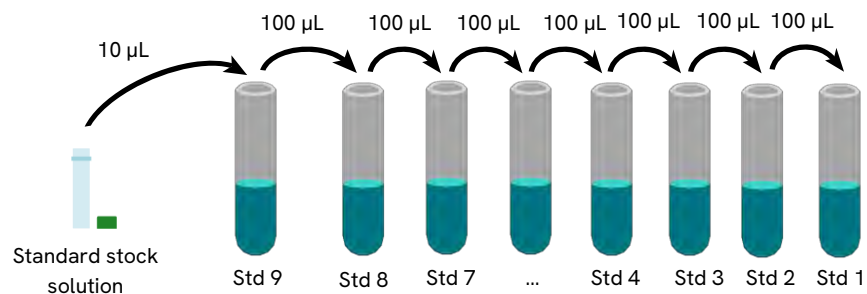
A recommended standard dilution procedure is listed and illustrated below.

- Dilute the standard stock solution 40-fold with diluent. This yields the high standard (Std 9 : 100 ng/mL) for the top of the curve. In practice:
  - e.g. take 10 $\mu\text{L}$  of the standard stock solution and add it to 390 $\mu\text{L}$  of diluent. Mix gently.
- Use the high standard (Std 9) to prepare the standard curve using 1/2 serial dilutions as follows:
  - Dispense 100 $\mu\text{L}$  of diluent in each vial from Std 8 to Std 1.
  - Add 100 $\mu\text{L}$  of standard 9 to 100 $\mu\text{L}$  of diluent, mix gently and repeat the 1/2 serial dilution to make standard solutions: 50, 25, 12.5, 6.25, 3.1, 1.6, 0.8, 0.4 ng/mL. This will create 9 standards for the analyte.
  - Std 0 (negative control) is diluent alone.

Step 1: dispense diluent in each vial.

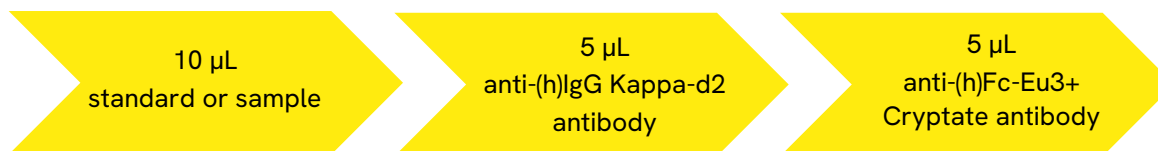


Step 2: dilute standards



## ASSAY PROTOCOL

Dispense the reagents in the following order:



The 2 HTRF® 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® compatible reader.

### For HTRF certified reader

For more information about HTRF® compatible readers and for set-up recommendations, please visit our website.

	Assays controls			
	Negative control	Cryptate control	Buffer control	Sample / Std
	<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	-	-	-	10 µL
Diluent	10 µL	10 µL	10 µL	-
Anti-(h) IgG Kappa-d2 antibody	5 µL	-	-	5 µL
Anti-(h) Fc-Eu <sup>3+</sup> -Cryptate antibody	5 µL	5 µL	-	5 µL
Detection buffer#3	-	5 µL	10 µL	-

## 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.

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

- 3) Calculate the % delta F which reflects the signal to background of the assay. The negative control plays the role of an internal assay control. Delta F is used for the comparison of day to day runs of the same assay.

$$\text{delta F (\%)} = \frac{\text{Ratio Standard or sample} - \text{Ratio Negative Control}}{\text{Ratio Negative Control}} \times 100$$

For more information about data reduction, please visit our website.

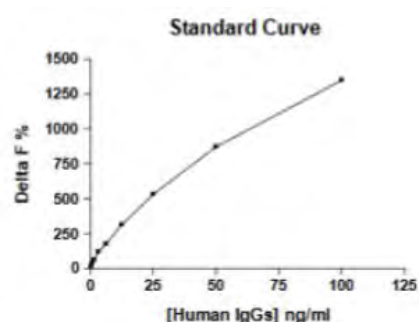
## RESULTS

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

Results may vary from one HTRF® 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)
Standard 0	Negative controle	948	2.0	0
Standard 1	0.4	1089	2.4	15
Standard 2	0.8	1298	2.8	37
Standard 3	1.6	1520	1.4	60
Standard 4	3.1	2066	2.9	118
Standard 5	6.25	2608	1.5	175
Standard 6	12.5	3921	1.0	314
Standard 7	25	5993	1.0	532
Standard 8	50	9220	1.2	872
Standard 9	100	13750	0.7	1350



## ASSAY CHARACTERISTICS

### Cross-reactivity

	Cross-reactivity %
Human Kappa	100
Human Lambda	0
Mouse Kappa	0
Human IgM (Kappa)	<1

### Detection limit

Human Kappa (IgG1) = 0.8 ng/mL

### 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 information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

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

revvity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.