

# MANUAL

Technology: HTRF<sup>®</sup> Cytokine

## HTRF High Performance Human IL-1 $\beta$ Detection Kit

Part number:	62HIL1B2PET
Test size	96 tests

Storage:  $\leq 60^{\circ}\text{C}$

Version: 1

Date: November 2023

## ASSAY PRINCIPLE

Revvity's human IL-1 $\beta$  assay is only intended for the quantitative measurement of IL-1 $\beta$  in supernatant using HTRF<sup>®</sup> technology. The assay is compatible with human samples, and is highly specific for IL-1 $\beta$ .

IL-1 $\beta$  is detected in a sandwich assay format using 2 different specific antibodies, one labeled with Europium Cryptate (donor) and the second with XL (acceptor).

The detection principle is based on HTRF<sup>®</sup> technology. When the labelled antibodies bind to the same antigen, 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). The two antibodies bind to the IL1 $\beta$  present in the sample, thereby generating FRET. Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the IL-1 $\beta$  concentration. (Fig. 1).

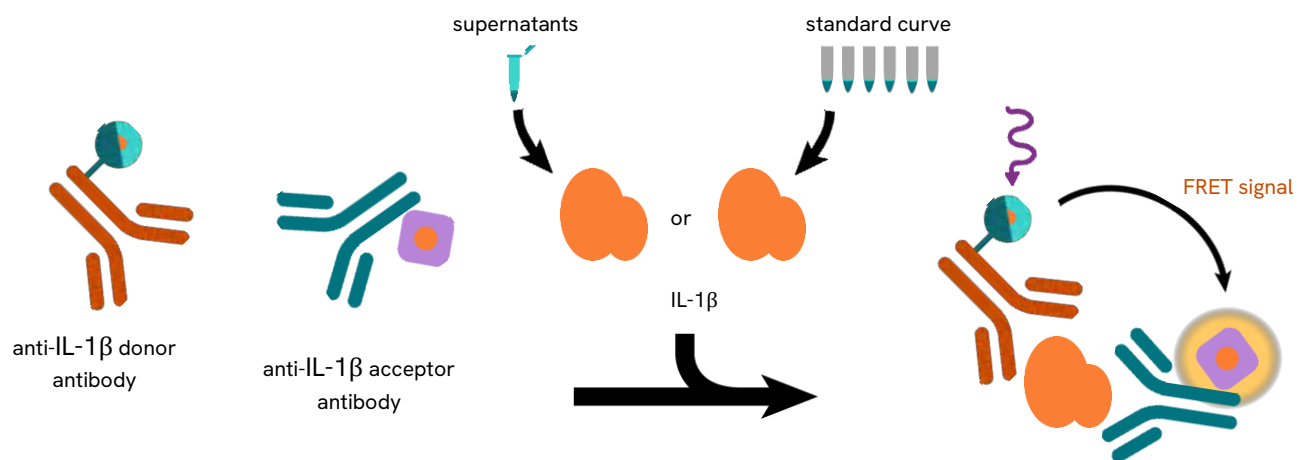
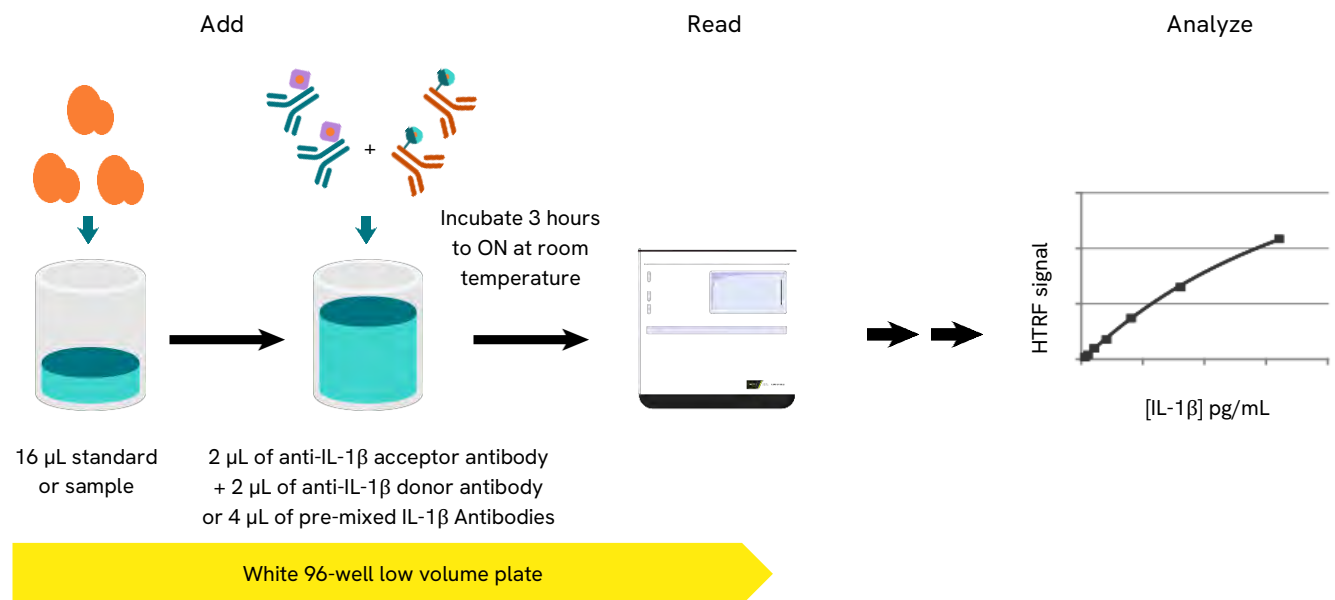


Figure 1: Principle of HTRF IL-1 $\beta$  sandwich assay

## PROTOCOL AT A GLANCE



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

## MATERIAL PROVIDED

KIT COMPONENTS	96 TESTS
IL-1 $\beta$ standard lyophilized	1 vial
IL-1 $\beta$ Eu cryptate antibody frozen 20X	1 vial 10 $\mu$ L
IL1 $\beta$ d2 antibody frozen 20X	1 vial 10 $\mu$ L
Diluent* #5 5X	1 vial 2 mL
Detection buffer** #3 Ready-to-use	1 vial 0.5 mL
Plate	1 plate HTRF 96-well low volume plate

\* To prepare working standard solutions, culture medium can be an alternative the diluent.

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

### Purchase separately

- HTRF®-Certified Reader. For a list of HTRF-compatible readers and set-up recommendations, please visit our website.
- 96-well or 384-well small volume (SV) detection microplates. For more information about microplate recommendations, please visit our website.

## 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.

Once reconstituted, standard stock solution may be frozen, and can be thawed only once.

Once thawed, antibody 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  $\leq -60^\circ\text{C}$ .

Volume of standard and antibody 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 the frozen reagents at room temperature.
- Before use, allow all kit's reagents to warm up at room temperature then
  - homogenize buffer and diluent with a vortex
  - centrifuge (NEVER vortex) the antibodies to gather all liquid at the bottom of the vial
- It is recommended to filter buffers before use.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.

**Take care to prepare stock and working solutions according to the directions for the kit size you have purchased.**

## To prepare reagent stock solutions

96 TESTS

### Anti-IL-1 $\beta$ Eu Cryptate antibody

Thaw the IL-1 $\beta$  Eu Cryptate antibody. Centrifuge.  
This stock solution can be frozen and stored at  $\leq -60^{\circ}\text{C}$



### Anti-IL-1 $\beta$ d2 antibody

Thaw the IL-1 $\beta$  XL antibody. Centrifuge.  
This stock solution can be frozen and stored at  $\leq -60^{\circ}\text{C}$ .



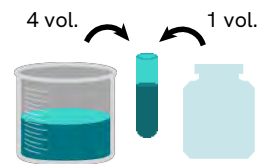
### IL-1 $\beta$ Standard

Reconstitute the IL-1 $\beta$  standard with distilled water.  
Volume of reconstitution is indicated on the vial label.  
The reconstituted standard solution can be frozen and stored at  $-60^{\circ}\text{C}$  or below.



### Diluent

Dilute 5-fold the 5 X diluent #5 with distilled water:  
homogenize the 5 X diluent #5 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.



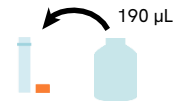
## To prepare antibody working solutions

Each well requires 4  $\mu\text{L}$  of pre-mixed IL-1 $\beta$  antibodies. Prepare the two antibody solutions in separate vials.

96 TESTS

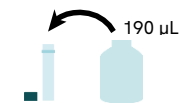
### IL-1 $\beta$ Eu Cryptate antibody

Dilute 20-fold the 20X stock solution (thawed reagent) of IL1 $\beta$  Eu Cryptate-antibody with detection buffer #3:  
Add 200 $\mu\text{L}$  of detection buffer directly in the thawed Eu Cryptate-antibody stock solution.



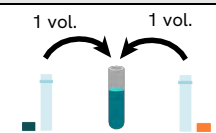
### IL-1 $\beta$ d2 antibody

Dilute 20-fold the 20X solution (thawed reagent) of IL1 $\beta$  d2 antibody with detection buffer #3:  
Add 200 $\mu\text{L}$  of detection buffer directly in thawed d2 antibody stock solution.



### Antibody Mix

Pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents: e.g. 200  $\mu\text{L}$  of d2 antibody + 200  $\mu\text{L}$  of Eu Cryptate antibody



## To prepare working standards solutions

- Each well requires 16  $\mu\text{L}$  of standard.
- Serially dilute the standard stock solution with with diluent #5. or with the cell culture medium used to prepare your samples supplemented with BSA or 10% FCS.
- **Due to the stability of the IL1 $\beta$ , it is mandatory to prepare the standard curve just before the assay.**
- 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 diluent.
- In order to counteract any standard sticking we recommend changing tips between each dilution.

## A recommended standard dilution procedure is listed and illustrated below:

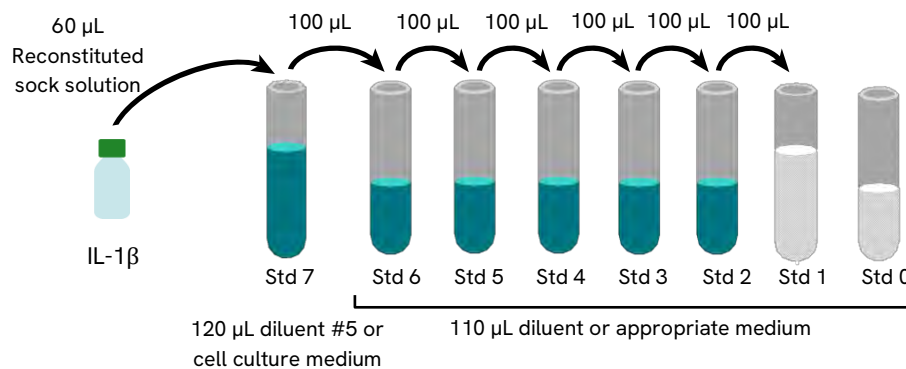
1. Reconstitute the standard vial with the volume indicated on the vial label using distilled water.
2. Prepare the following dilutions:

- Dilute the reconstituted standard stock solution 3-fold with diluent or with cell culture medium.

In practice: take 60  $\mu\text{L}$  of stock solution and add it to 120  $\mu\text{L}$  of diluent or cell culture medium. Mix gently. This yields the high standard (Std 7: 6500 pg/mL) for the top of the curve.

- Use the high standard (Std 7) to prepare the standard curve using serial dilutions as follows:
  - Dispense 110  $\mu\text{L}$  of diluent or cell culture medium into each vial from Std 6 to Std 0
  - Add 100  $\mu\text{L}$  of standard to 110  $\mu\text{L}$  of diluent or cell culture medium, mix gently and repeat the 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 diluent or appropriate culture medium alone








STANDARD	SERIAL DILUTIONS	WORKING SOLUTIONS
Standard Stock solution	Reconstitute the vial following the indications given on the vial label	19.5 ng/mL
Standard 7	60 $\mu\text{L}$ Standard stock Solution + 120 $\mu\text{L}$ diluent	6 500 pg/mL
Standard 6	100 $\mu\text{L}$ standard 7 + 110 $\mu\text{L}$ diluent	3095.2 pg/mL
Standard 5	100 $\mu\text{L}$ standard 6 + 110 $\mu\text{L}$ diluent	1473.9 pg/mL
Standard 4	100 $\mu\text{L}$ standard 5 + 110 $\mu\text{L}$ diluent	701.9 pg/mL
Standard 3	100 $\mu\text{L}$ standard 4 + 110 $\mu\text{L}$ diluent	334.2 pg/mL
Standard 2	100 $\mu\text{L}$ standard 3 + 110 $\mu\text{L}$ diluent	159.1 pg/mL
Standard 1	100 $\mu\text{L}$ standard 2 + 110 $\mu\text{L}$ diluent	75.8 pg/mL
Standard 0	110 $\mu\text{L}$ diluent	0

## To prepare samples

- Each well requires 16  $\mu\text{L}$  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 (Std 7) must be diluted diluent #5 in your appropriate sample medium.

## ASSAY PROTOCOL

		STANDARD (STD 0 – STD 7)	SAMPLES
<b>Step 1</b>		Dispense 16 $\mu\text{L}$ of each IL-1 $\beta$ standard (Std 0 - Std 7) into each standard well	Dispense 16 $\mu\text{L}$ of each sample into each sample well
<b>Step 2</b>		Add 2 $\mu\text{L}$ of IL-1 $\beta$ d2 antibody working solution to all wells	
<b>Step 3</b>		Add 2 $\mu\text{L}$ of IL-1 $\beta$ Eu Cryptate antibody working solution to all wells.	
<b>Step 4</b>		Seal the plate and incubate from 3 hours at RT	
<b>Step 5</b>		Remove the plate sealer and read on an HTRF® compatible reader	

	1	2	3	4	5	6
A	16 $\mu\text{L}$ Std 0 (Negative control) 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well A1	Repeat Well A1	16 $\mu\text{L}$ sample 1 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well A4	Repeat Well A4
B	16 $\mu\text{L}$ Std 1 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well B1	Repeat Well B1	16 $\mu\text{L}$ sample 2 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well B4	Repeat Well B4
C	16 $\mu\text{L}$ Std 2 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well C1	Repeat Well C1	16 $\mu\text{L}$ sample 3 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well C4	Repeat Well C4
D	16 $\mu\text{L}$ Std 3 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well D1	Repeat Well D1	16 $\mu\text{L}$ sample... 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well D4	Repeat Well D4
E	16 $\mu\text{L}$ Std 4 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well E1	Repeat Well E1	16 $\mu\text{L}$ sample... 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well E4	Repeat Well E4
F	16 $\mu\text{L}$ Std 5 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well F1	Repeat Well F1	16 $\mu\text{L}$ sample... 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well F4	Repeat Well F4
G	16 $\mu\text{L}$ Std 6 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well G1	Repeat Well G1	16 $\mu\text{L}$ sample... 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well G4	Repeat Well G4
H	16 $\mu\text{L}$ Std 7 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well H1	Repeat Well H1	16 $\mu\text{L}$ sample... 2 $\mu\text{L}$ IL1 $\beta$ -d2 2 $\mu\text{L}$ IL1 $\beta$ -Eu Cryptate	Repeat Well H4	Repeat Well H4

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A																								
B																								
C																								
D																								
E																								
F																								
G																								
H																								
I																								
J																								
K																								
L																								
M																								

## DATA REDUCTION & INTERPRETATION

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

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

- 2) Calculate the delta ratio of the acceptor and donor emission signals for each individual well. The Standard 0 (Negative control) plays the role of an internal assay control.

$$\text{Delta Ratio} = \text{Ratio Standard or sample} - \text{Ratio Standard 0}$$

- 3) 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 our website.

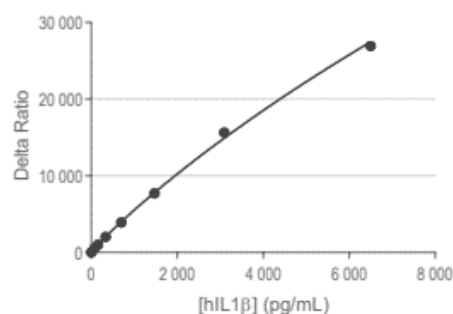
## RESULTS

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 with  $1/Y^2$ ) model

		Ratio (1)	CV% (2)
Standard 0	Negative control	1127	7.5%
Standard 1	75.8 pg/mL	1942	1.8%
Standard 2	159.1 pg/mL	2905	3.8%
Standard 3	334.2 pg/mL	4526	0.8%
Standard 4	701.9 pg/mL	7930	1.5%
Standard 5	1 473.9 pg/mL	14276	4.9%
Standard 6	3 095.2 pg/mL	22537	2.4%
Standard 7	6 500 pg/mL	29746	2.4%



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.



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.