

MANUAL

Technology: HTRF™ Cytokine

HTRF High Performance Human IL-6 Kit

Part number:	62HIL6V2PEG	62HIL6V2PEH
Test size	500 tests	10,000 tests

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

Version: 01

Date: July 2024

ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of human IL-6 in supernatant and offers a fast alternative to ELISA.

The detection principle of this kit is based on HTRF™ technology (Homogeneous Time-Resolved Fluorescence). As shown in Figure 1, IL-6 is detected in a sandwich assay by using anti-IL-6 antibody labeled with Europium cryptate (donor), and anti-IL-6 antibody labeled 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 (665 nm). Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the IL-6 concentration.

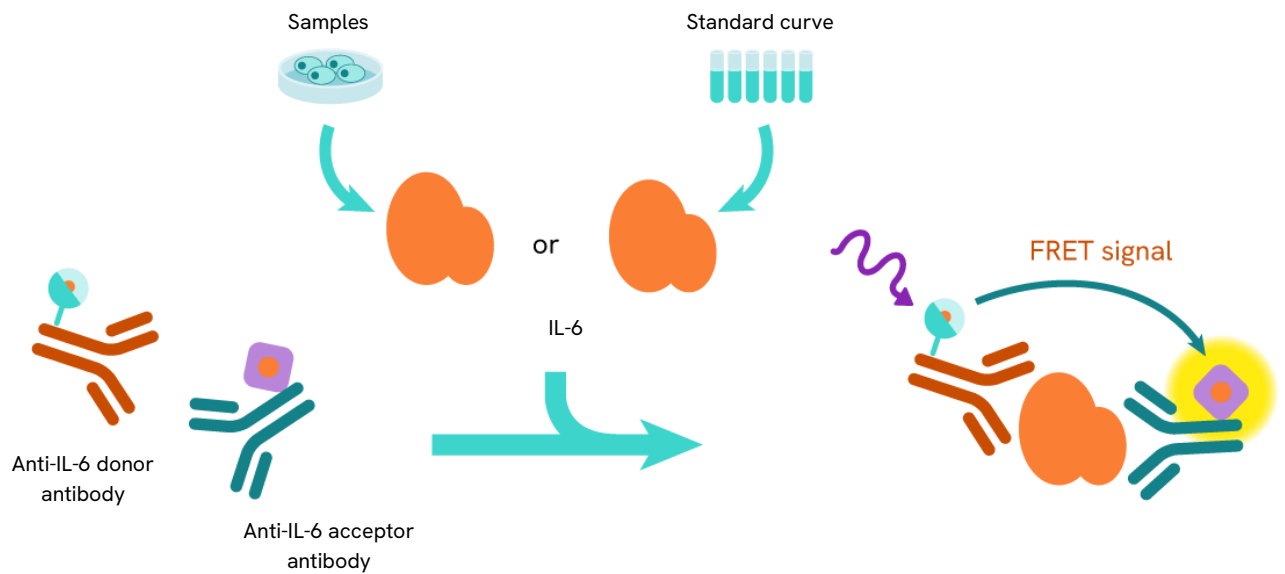
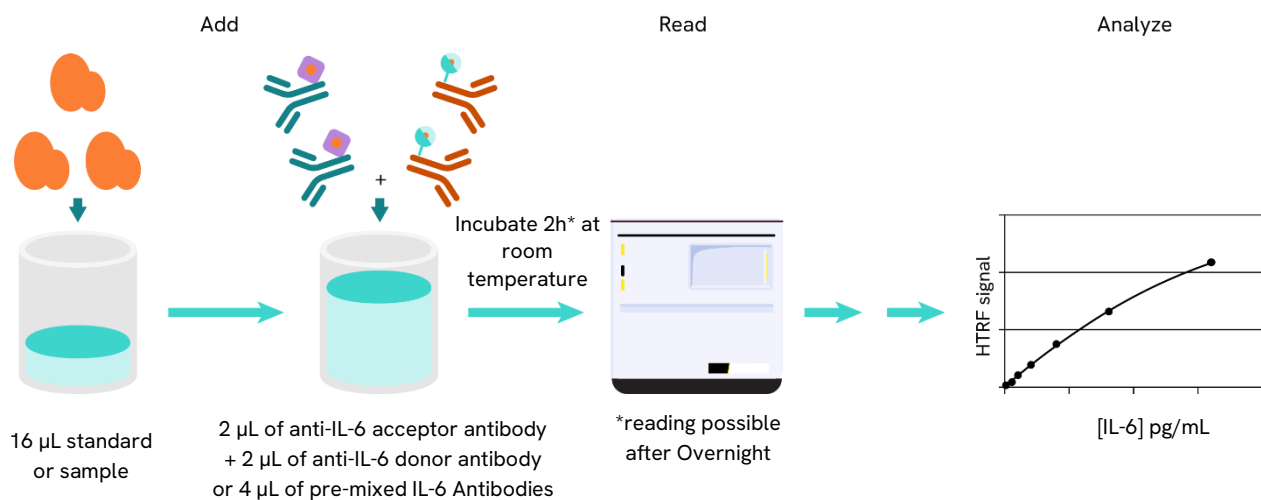



Figure 1: Principle of HTRF IL-6 sandwich assay

PROTOCOL AT A GLANCE



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

MATERIAL PROVIDED

KIT COMPONENTS	500 TESTS*			10,000 TESTS*		
IL-6 Standard Lyophilized		green cap	2 vials		green cap	2 vials
IL-6 Eu Cryptate Antibody Frozen 20X		orange cap	1 vial 50 µL		red cap	1 vial 1 mL
IL-6 d2 Antibody Frozen 20X		blue cap	1 vial 50 µL		purple cap	1 vial 1 mL
Diluent** #5 5X		white cap	1 vial 2 mL		white cap	1 vial 10 mL
Detection Buffer*** #3 Ready-to-use		red cap	2 vials 1.5 mL		red cap	1 vial 50 mL

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

** Medium like cell culture medium can be an alternative to the diluent.

*** The Detection buffer is used to prepare working solutions of acceptor and donor reagents.

Purchase separately

- HTRF™-Certified Reader. **Make sure the setup for Eu Cryptate is used.** For a list of HTRF-compatible readers and set-up recommendations, please visit our website.
- Small volume (SV) detection microplates. Use white plate only. For more information about microplate recommendations, please visit our website.

STORAGE AND STABILITY

Kit

- Store the kit at -20°C or below.
- Under proper storage conditions, reagents are stable until the expiry date indicated on the label.
- Diluent and detection buffer are shipped frozen but can be stored at 2-8°C in your premises.

Reagents

Once reconstituted, standard stock solution may be frozen at ≤-60°C, 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 ≤-60°C for standard and at ≤-20°C for antibodies.

Volume of standard and antibody aliquots should not be under 10 µL.

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, allow them to warm up to room temperature for at least 30 mins before use.
- Before use, allow Diluent and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- IL-6 standards (for standard curve) must be prepared in diluent or in the same medium as the samples.

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

To prepare reagent stock solutions

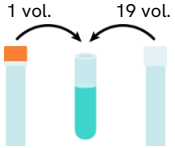
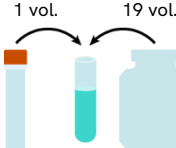
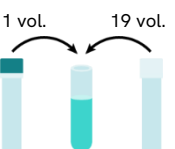

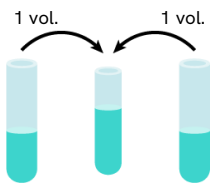
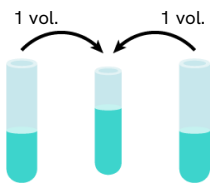
500 TESTS		10,000 TESTS	
Anti-IL-6 Eu Cryptate antibody			
Thaw the IL-6 Eu Cryptate antibody. Mix gently. This 20X stock solution can be frozen and stored at -20°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -20°C or below.			Thaw the IL-6 Eu Cryptate antibody. Mix gently. This 20X stock solution can be frozen and stored at -20°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -20°C or below.
Anti-IL-6 d2 antibody			
Thaw the IL-6 d2 antibody. Mix gently. This 20X stock solution can be frozen and stored at -20°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -20°C or below.			Thaw the IL-6 d2 antibody. Mix gently. This 20X stock solution can be frozen and stored at -20°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -20°C or below.
IL-6 Standard			
Reconstitute the IL-6 standard with distilled water. Volume of reconstitution is indicated on the vial label. The reconstituted standard solution can be frozen and stored at -60°C or below.			Reconstitute the IL-6 standard with distilled water. Volume of reconstitution is indicated on the vial label. The reconstituted standard solution can be frozen and stored at -60°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. This 1X diluent can be frozen and stored at -20°C or below.			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., 10 mL of diluent + 40 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -20°C or below.
Detection buffer			

The Detection buffer is ready-to-use.

To prepare antibody working solutions

Each well requires 2 μL of IL-6-Eu Cryptate Antibody and 2 μL of IL-6 d2 Antibody.

Prepare the two antibody solutions in separate vials.

500 TESTS		10,000 TESTS	
IL-6 Eu Cryptate antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of IL-6 Eu Cryptate antibody with Detection buffer #3: add 1 volume of Eu Cryptate antibody stock solution in 19 volumes of detection buffer (e.g. 10 μL of Eu Cryptate antibody stock solution + 190 μL of detection buffer).			Dilute 20-fold the 20X stock solution (thawed reagent) of IL-6 Eu Cryptate antibody with Detection buffer #3: add 1 volume of Eu Cryptate antibody stock solution in 19 volumes of detection buffer (e.g. 10 μL of Eu Cryptate antibody stock solution + 190 μL of detection buffer).
IL-6 d2 antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of IL-6 d2 antibody with Detection buffer #3: add 1 volume of d2 antibody stock solution in 19 volumes of detection buffer (e.g. 10 μL of Eu Cryptate antibody stock solution + 190 μL of detection buffer).			Dilute 20-fold the 20X stock solution (thawed reagent) of IL-6 d2 antibody with Detection buffer #3: add 1 volume of d2 antibody stock solution in 19 volumes of detection buffer (e.g. 10 μL of Eu Cryptate antibody stock solution + 190 μL of detection buffer).
Antibody Mix			
It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).			It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 20 mL of d2 antibody + 20 mL of Cryptate antibody).

To prepare working standards solutions

- Each well requires 16 μL of standard.
- Dilute the standard stock solution serially with diluent #5 (1X)
- If culture medium is used to dilute the standard, we recommend to supplement it with serum (2 to 10%) or BSA (0.2 to 1%) in order to avoid IL-6 sticking to assay plates.
- 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 #5 (1X).
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

A recommended standard dilution procedure is listed and illustrated below:

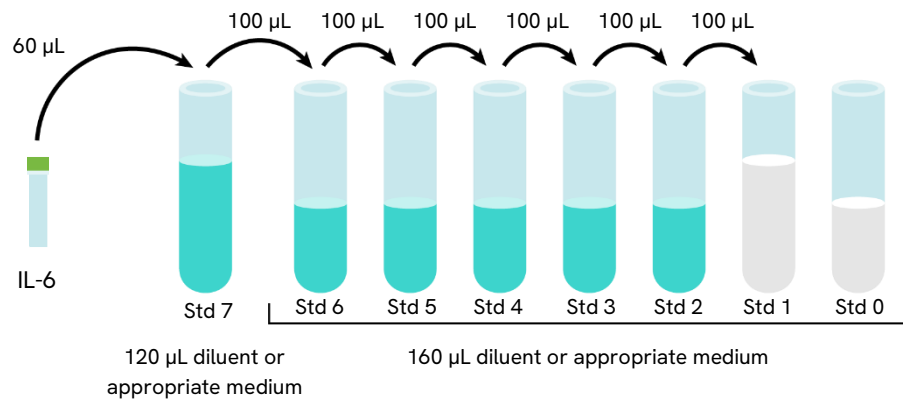
Dilute the standard stock solution 3-fold with diluent; this yields the Standard Max solution (12 000 pg/mL)

Dilute the standard stock solution 3-fold with diluent #5 (1X) to prepare high standard (Std 7): e.g. take 60 µL of standard stock solution and add it to 120 µL of diluent #5 (1X). Mix gently.

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

- Dispense 160 µL of diluent #5 (1X) in each vial from Std 6 to Std 0.
- Add 100 µL of standard to 160 µL of diluent #5 (1X), 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 #5 (1X) or appropriate culture medium alone








STANDARD	SERIAL DILUTIONS	WORKING SOLUTIONS
Standard Stock solution	Reconstituted lyophilizate	36 ng/mL
Standard 7	60 µL Standard stock Solution + 120 µL diluent 1X	12 000 pg/mL
Standard 6	100 µL standard 7 + 160 µL diluent 1X	4615.4 pg/mL
Standard 5	100 µL standard 6 + 160 µL diluent 1X	1775.1 pg/mL
Standard 4	100 µL standard 5 + 160 µL diluent 1X	682.7 pg/mL
Standard 3	100 µL standard 4 + 160 µL diluent 1X	262.6 pg/mL
Standard 2	100 µL standard 3 + 160 µL diluent 1X	101 pg/mL
Standard 1	100 µL standard 2 + 160 µL diluent 1X	38.8 pg/mL
Standard 0	160 µL diluent 1X	0

To prepare samples

- Each well requires 16 µL 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 in your appropriate sample medium.
- To obtain additional information or support, please contact the HTRF technical support team.

ASSAY PROTOCOL

		STANDARD (STD 0 – STD 7)	SAMPLES
Step 1		Dispense 16 µL of each IL-6 standard (Std 0 - Std 7) into each standard well	Dispense 16 µL of each sample into each sample well
Step 2		Add 2 µL of IL-6 d2 antibody working solution to all wells	
Step 3		Add 2 µL of IL-6 Eu Cryptate antibody working solution to all wells.	
Step 4		Seal the plate and incubate from 2 hours at RT *Reading possible after Overnight	
Step 5		Remove the plate sealer and read on an HTRF™ compatible reader	

	1	2	3	4	5	6
A	16 µL Std 0 (Negative control) 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well A1	Repeat Well A1	16 µL sample 1 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well A4	Repeat Well A4
B	16 µL Std 1 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well B1	Repeat Well B1	16 µL sample 2 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well B4	Repeat Well B4
C	16 µL Std 2 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well C1	Repeat Well C1	16 µL sample 3 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well C4	Repeat Well C4
D	16 µL Std 3 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well D1	Repeat Well D1	16 µL sample... 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well D4	Repeat Well D4
E	16 µL Std 4 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well E1	Repeat Well E1	16 µL sample... 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well E4	Repeat Well E4
F	16 µL Std 5 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well F1	Repeat Well F1	16 µL sample... 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well F4	Repeat Well F4
G	16 µL Std 6 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well G1	Repeat Well G1	16 µL sample... 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well G4	Repeat Well G4
H	16 µL Std 7 2 µL IL-6-d2 2 µL IL-6-Eu Cryptate	Repeat Well H1	Repeat Well H1	16 µL sample... 2 µL IL-6-d2 2 µL IL-6-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																								
N																								
O																								
P																								
Q																								

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 our website.

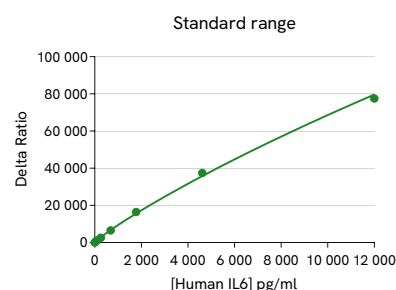
RESULTS

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

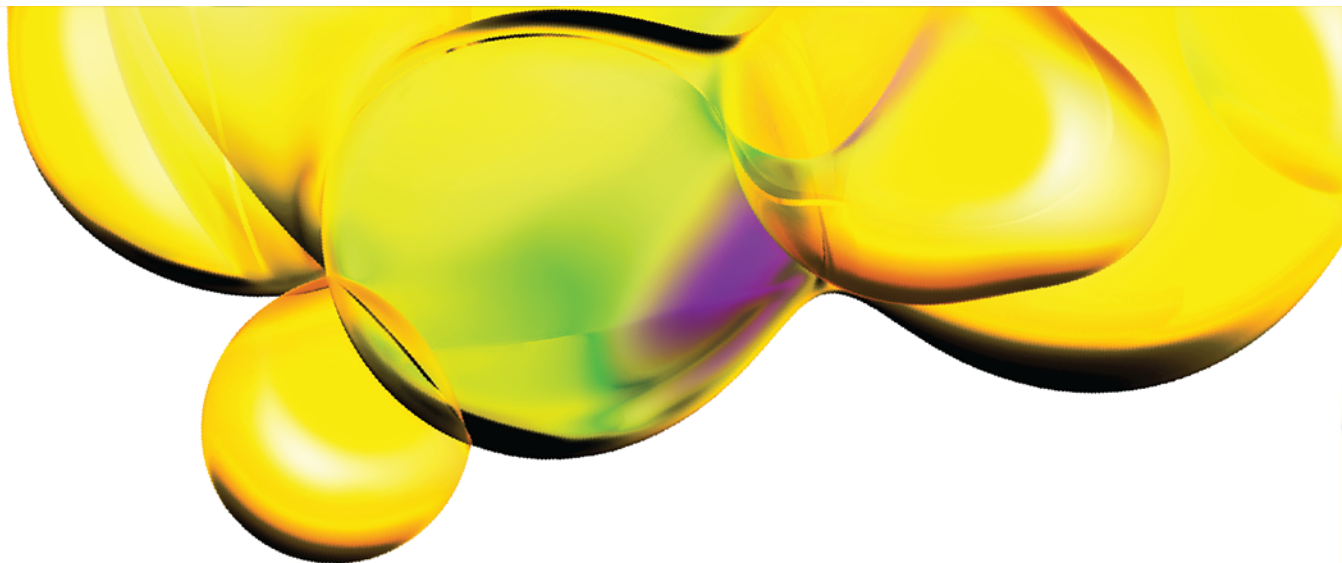
		Ratio (1)	CV% (2)
Standard 0	Negative control	1117	8%
Standard 1	38.8 pg/mL	1777	5,1%
Standard 2	101 pg/mL	2409	4,4%
Standard 3	262.6 pg/mL	3831	0,5%
Standard 4	682.7 pg/mL	7714	3,2%
Standard 5	1775.1 pg/mL	17712	2,1%
Standard 6	4615.4 pg/mL	38652	1,2%
Standard 7	12 000 pg/mL	78740	2,6%



REACH European regulations and compliance

This product and/or some of its components may 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

revvity

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

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.