

# MANUAL

Technology: HTRF™ Cytokine

# HTRF High Performance Human TNF alpha Kit

|              |              |
|--------------|--------------|
| Part number: | 62HTNFAV2PET |
| Test size    | 96 tests     |

Storage:  $\leq -16^{\circ}\text{C}$

Version: 01

Date: January 2025

## ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of human TNF $\alpha$  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, TNF $\alpha$  is detected in a sandwich assay by using anti- TNF $\alpha$  antibody labeled with Europium cryptate (donor), and anti- TNF $\alpha$  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 TNF $\alpha$  concentration.

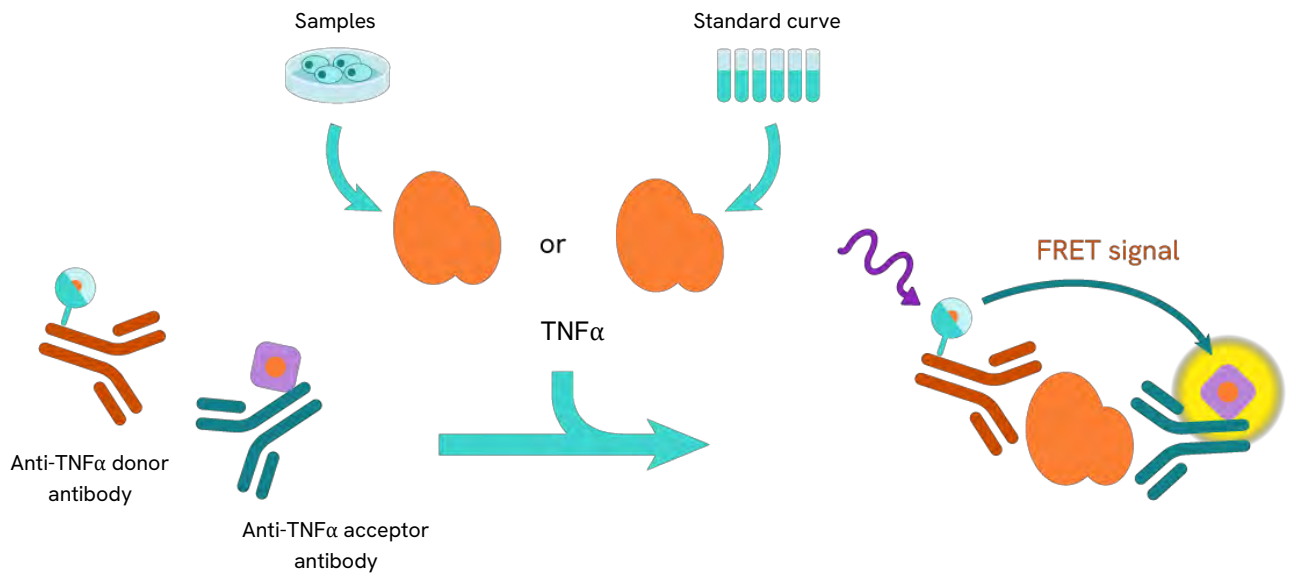
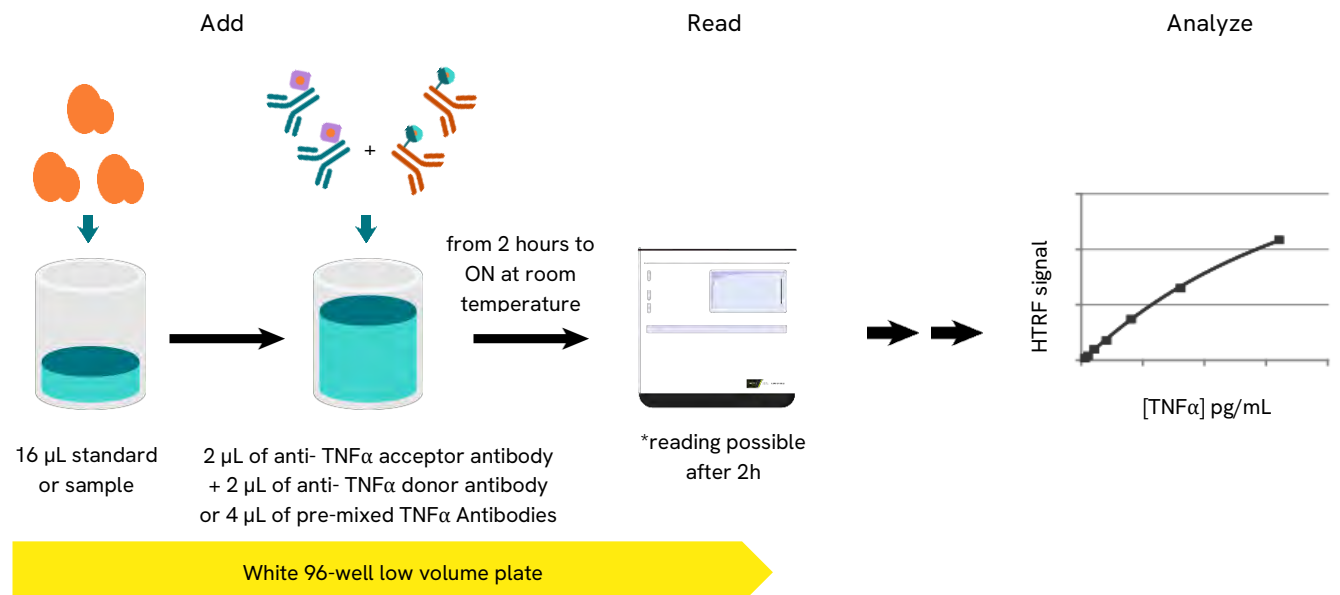







Figure 1: Principle of HTRF TNF $\alpha$  sandwich assay

## PROTOCOL AT A GLANCE



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

## MATERIAL PROVIDED

| KIT COMPONENTS                               | 96 TESTS  |                 |  |
|--|---|-----------------|--|
| TNF $\alpha$ Standard Lyophilized            |  | green cap       | 1 vial                                   |
| TNF $\alpha$ Eu Cryptate Antibody Frozen 20X |  | orange cap      | 1 vial<br>10 $\mu$ L                     |
| TNF $\alpha$ d2 Antibody Frozen 20X          |  | blue cap        | 1 vial<br>10 $\mu$ L                     |
| Diluent* #1                                  |  | Transparent cap | 1 vial<br>20 mL                          |
| Detection Buffer** #3 Ready-to-use           |  | Transparent cap | 1 vial<br>0.5 mL                         |
| Plate  |   |                 | 1 plate<br>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 -16°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  $\leq -16^{\circ}\text{C}$ .

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.



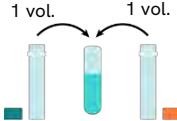
**Take care to prepare stock and working solutions according to the directions for the kit size you have purchased.**

## To prepare reagent stock solutions

| 96 TESTS  |   |
|---|---|
| <b>Anti-TNF<math>\alpha</math> Eu Cryptate antibody</b>   |   |
| Thaw the TNF $\alpha$ Eu Cryptate antibody. Centrifuge.<br>This stock solution can be frozen and stored at $\leq -16^{\circ}\text{C}$   |  |
| <b>Anti-TNF<math>\alpha</math> d2 antibody</b>  |   |
| Thaw the TNF $\alpha$ d2 antibody. Centrifuge.<br>This stock solution can be frozen and stored at $\leq -16^{\circ}\text{C}$ .  |  |
| <b>TNF<math>\alpha</math> Standard</b>  |   |
| Reconstitute the TNF $\alpha$ standard with distilled water.<br>Volume of reconstitution is indicated on the vial label.<br>The reconstituted standard solution can be frozen and stored at $-60^{\circ}\text{C}$ or below. |  |
| <b>Diluent</b>  |   |
| The Diluent is ready-to-use.  |   |

## To prepare antibody working solutions

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

| 96 TESTS   |   |
|--|---|
| <b>TNF<math>\alpha</math> Eu Cryptate antibody</b>   |   |
| Dilute the stock solution (thawed reagent) of TNF $\alpha$ Eu Cryptate-antibody with detection buffer #3:<br>Add 200 $\mu\text{L}$ of detection buffer directly in the thawed Eu Cryptate-antibody stock solution. |  |
| <b>TNF<math>\alpha</math> d2 antibody</b>  |   |
| Dilute the stock solution (thawed reagent) of TNF $\alpha$ d2 antibody with detection buffer #3:<br>Add 200 $\mu\text{L}$ of detection buffer directly in thawed d2 antibody stock solution.                       |  |
| <b>Antibody Mix</b>  |   |
| Pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents: e.g. 210 $\mu\text{L}$ of d2 antibody + 210 $\mu\text{L}$ 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 diluent #1 or with the cell culture medium used to prepare your samples supplemented with BSA or 10% FCS.
- **Due to the stability of the  $\text{TNF}\alpha$ , 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:

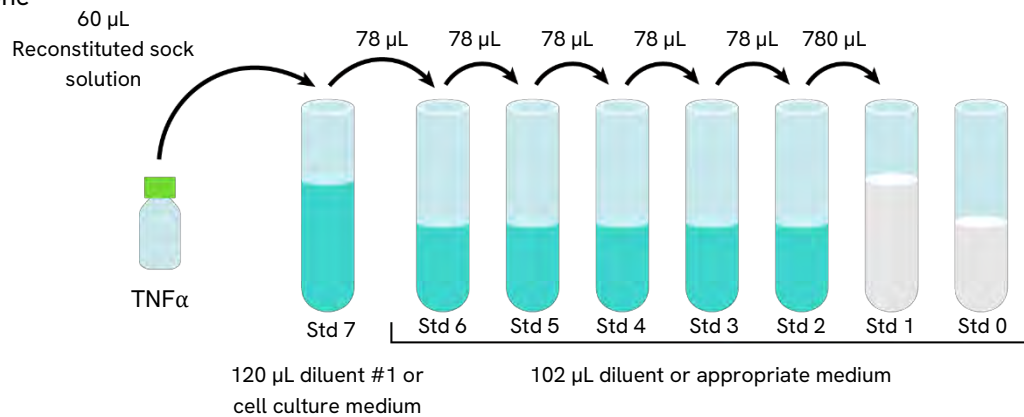
1. Reconstitute the standard vial with the volume indicated on the vial label using distilled water.
2. Prepare the following dilutions:

- 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: 5 500  $\text{pg}/\text{mL}$ ) for the top of the curve.

- Use the high standard (Std 7) to prepare the standard curve using serial dilutions as follows:
  - Dispense 102  $\mu\text{L}$  of diluent or cell culture medium into each vial from Std 6 to Std 0
  - Add 78  $\mu\text{L}$  of standard to 102  $\mu\text{L}$  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 | Reconstituted lyophilisate  | 16.5 $\text{ng}/\text{mL}$  |
| Standard 7              | 60 $\mu\text{L}$ Standard stock Solution + 120 $\mu\text{L}$ diluent #1 | 5 500 $\text{pg}/\text{mL}$ |
| Standard 6              | 78 $\mu\text{L}$ standard 7 + 102 $\mu\text{L}$ diluent #1              | 2391 $\text{pg}/\text{mL}$  |
| Standard 5              | 78 $\mu\text{L}$ standard 6 + 102 $\mu\text{L}$ diluent #1              | 1040 $\text{pg}/\text{mL}$  |
| Standard 4              | 78 $\mu\text{L}$ standard 5 + 102 $\mu\text{L}$ diluent #1              | 452 $\text{pg}/\text{mL}$   |
| Standard 3              | 78 $\mu\text{L}$ standard 4 + 102 $\mu\text{L}$ diluent #1              | 196 $\text{pg}/\text{mL}$   |
| Standard 2              | 78 $\mu\text{L}$ standard 3 + 102 $\mu\text{L}$ diluent #1              | 85 $\text{pg}/\text{mL}$    |
| Standard 1              | 78 $\mu\text{L}$ standard 2 + 102 $\mu\text{L}$ diluent #1              | 37 $\text{pg}/\text{mL}$    |
| Standard 0              | 102 $\mu\text{L}$ diluent #1  | 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 #1 or in your appropriate sample medium.

## ASSAY PROTOCOL

|        |   | STANDARD (STD 0 – STD 7)  | SAMPLES  |
|--------|---|---|--|
| Step 1 |  | Dispense 16 $\mu\text{L}$ of each TNF $\alpha$ standard (Std 0 - Std 7) into each standard well | Dispense 16 $\mu\text{L}$ of each sample into each sample well |
| Step 2 |  | Add 2 $\mu\text{L}$ of TNF $\alpha$ d2 antibody working solution to all wells                   |  |
| Step 3 |  | Add 2 $\mu\text{L}$ of TNF $\alpha$ Eu Cryptate antibody working solution to all wells.         |  |
| Step 4 |  | Seal the plate and incubate at RT<br>*Reading possible after 2h                                 |  |
| Step 5 |  | 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)<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate | Repeat Well A1 | Repeat Well A1 | 16 $\mu\text{L}$ sample 1<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate  | Repeat Well A4 | Repeat Well A4 |
| B | 16 $\mu\text{L}$ Std 1<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well B1 | Repeat Well B1 | 16 $\mu\text{L}$ sample 2<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate  | Repeat Well B4 | Repeat Well B4 |
| C | 16 $\mu\text{L}$ Std 2<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well C1 | Repeat Well C1 | 16 $\mu\text{L}$ sample 3<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate  | Repeat Well C4 | Repeat Well C4 |
| D | 16 $\mu\text{L}$ Std 3<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well D1 | Repeat Well D1 | 16 $\mu\text{L}$ sample...<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate | Repeat Well D4 | Repeat Well D4 |
| E | 16 $\mu\text{L}$ Std 4<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well E1 | Repeat Well E1 | 16 $\mu\text{L}$ sample...<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate | Repeat Well E4 | Repeat Well E4 |
| F | 16 $\mu\text{L}$ Std 5<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well F1 | Repeat Well F1 | 16 $\mu\text{L}$ sample...<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate | Repeat Well F4 | Repeat Well F4 |
| G | 16 $\mu\text{L}$ Std 6<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well G1 | Repeat Well G1 | 16 $\mu\text{L}$ sample...<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate | Repeat Well G4 | Repeat Well G4 |
| H | 16 $\mu\text{L}$ Std 7<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-Eu Cryptate                    | Repeat Well H1 | Repeat Well H1 | 16 $\mu\text{L}$ sample...<br>2 $\mu\text{L}$ TNFa-d2<br>2 $\mu\text{L}$ TNFa-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 |   |   |   |   |   |   |   |   |   |    |    |    |    |    |    |    |    |    |    |    |    |    |    |    |

## 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 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.

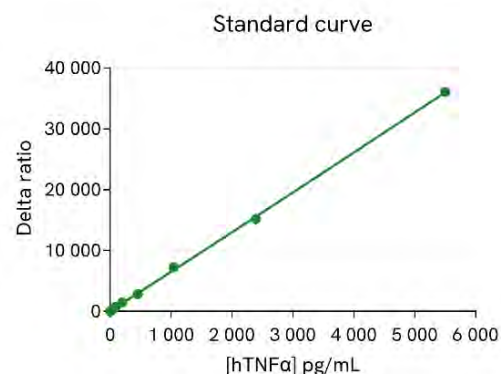
## RESULTS

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) |
|-------------------|-------------------------|-----------|---------|
| <b>Standard 0</b> | <b>Negative control</b> | 1171      | 3,1%    |
| Standard 1        | 37 pg/mL                | 1407      | 3,4%    |
| Standard 2        | 85 pg/mL                | 1878      | 8,0%    |
| Standard 3        | 196 pg/mL               | 2610      | 6,7%    |
| Standard 4        | 452 pg/mL               | 4273      | 3,6%    |
| Standard 5        | 1 040 pg/mL             | 8438      | 2,6%    |
| Standard 6        | 2391 pg/mL              | 16353     | 2,9%    |
| Standard 7        | 5 500 pg/mL             | 37235     | 0,8%    |



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