

HTRF® Research Reagents

### HTRF Human Granzyme B kit - 500 pts

Product Number : 62HGRZBPEG Lot Number : 07H

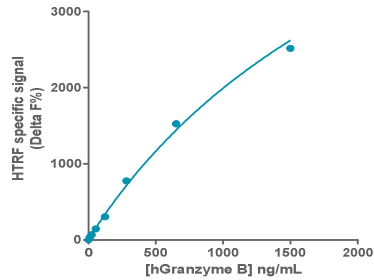
Expiration date 16/08/2024 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	29A
HTRF Human Granzyme B kit - Std	04A
HTRF Human Granzyme B kit - 500 pts d2 antibody	07A
HTRF Human Granzyme B kit - 500 pts Eu Cryptate antibody	07A
HTRF Human Granzyme B kit - modop	5

**Results :**

Incubation : **Overnight at RT**  
 Plate : **384 wells small volume white plate 20 µL**  
 Reader : **Pherastar Plus**  
 Fitting model : **4 Parameter Logistic (4PL)**

	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
<b>Std 0</b>	0	0%	0-0
<b>Std 1</b>	10	32	15-48
<b>Std 2</b>	23	68	60-93
<b>Std 3</b>	54	147	80-237
<b>Std 4</b>	123	305	321-414
<b>Std 5</b>	283	778	668-1112
<b>Std 6</b>	852	1527	1295-2145
<b>Std 7</b>	1500	2519	2194-3101



**Assay characteristics :**

	Found value	Specifications Rev 01
	/	/
	/	/
	/	/
	/	/

HTRF Human Granzyme B kit - 500 pts - Lot n° : 07H

Accepted

Rejected

Quality control date : 2023-09-01	Supervisor's initials : PL
-----------------------------------	----------------------------

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.