

HTRF® Research Reagents

HTRF Human Granzyme B kit - 500 pts

Product Number : 62HGRZBPEG Lot Number : 07K

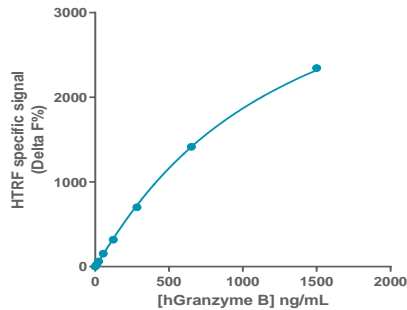
Expiration date : 2024-08-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	19B
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human Granzyme B kit - Std	04RA
HTRF Human Granzyme B kit - 500 pts d2 antibody	06BR
HTRF Human Granzyme B kit - 500 pts Eu Cryptate antibody	07RA
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
Std 0	0	0%	0-0
Std 1	10	23	15-48
Std 2	23	62	60-93
Std 3	54	154	80-237
Std 4	123	319	321-414
Std 5	283	701	668-1112
Std 6	852	1416	1295-2145
Std 7	1500	2345	2194-3101



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

Accepted Rejected

Quality control date : 2023-11-13	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.