

HTRF® Research Reagents

HTRF Human Granzyme B kit - 500 pts

Product Number : 62HGRZBPEG Lot Number : 07M

Expiration date : 2026-08-16 Document version : 1

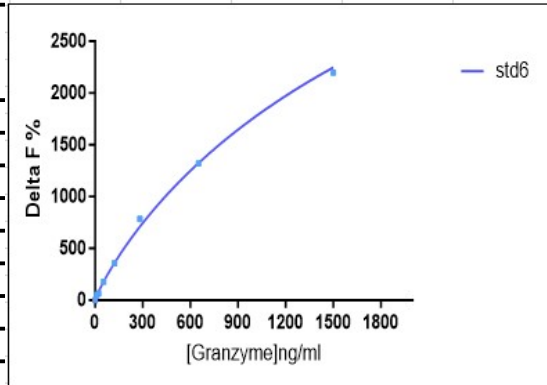
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human Granzyme B kit - Std	06A
HTRF Human Granzyme B kit - 500 pts d2 antibody	06BR
HTRF Human Granzyme B kit - 500 pts Eu Cryptate antibody	07RA

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	10	48	15-48
Std 2	23	66	60-93
Std 3	54	179	80-237
Std 4	123	358	321-414
Std 5	283	784	668-1112
Std 6	852	1319	1295-2145
Std 7	1500	2197	2194-3101



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

Accepted

Rejected

Quality control date : 2024-04-16	Supervisor's initials : AC
-----------------------------------	----------------------------

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.