

HTRF® Research Reagents

HTRF Human Progranulin kit - 500 pts

Product Number : 62PROPEG Lot Number : 02A

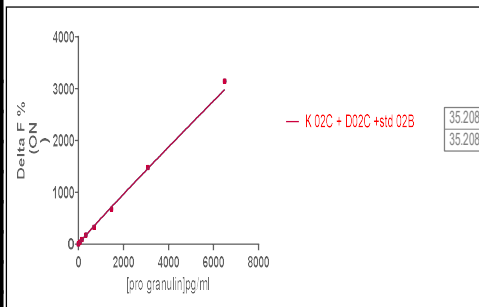
Expiration date : 2026-06-07 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21RA
HTRF Detect. Buf. 3 - 1.5 mL		30AR
HTRF Human Progranulin kit - Std		02B
HTRF Human Progranulin kit - 500 pts d2 antibody		02C
HTRF Human Progranulin kit - 500 pts Eu Cryptate antibody		02C

Results :

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

hProgranulin	Conc (pg/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	338	42%	39-67
Std 2	777	87%	83-137
Std 3	1787	178%	163-258
Std 4	4109	325%	316-471
Std 5	9452	677%	612-914
Std 6	21739	1481%	1286-1757
Std 7	50000	3145%	2358-3089



Assay characteristics :

	Found value	Specifications Rev 01
St 7 / St 6	2.12	>1.6

HTRF Human Progranulin kit - 500 pts - Lot n° : 02A

Accepted

Rejected

Quality control date :	2024-05-23	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.