

HTRF® Research Reagents

HTRF (h/m) HTT Mutant Kit - 500 pts

Product Number : 64HTTMPEG Lot Number : 03C

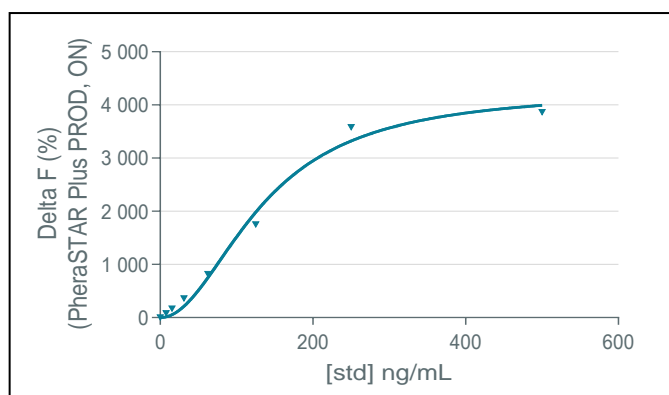
Expiration date : 2026-06-07 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 15 - 2 mL	09A
HTRF (h/m) HTT Mutant Kit - 500 pts d2 antibody	03A
HTRF (h/m) HTT Mutant Kit - 500 pts Tb Cryptate antibody	03A
HTRF P-T prot. - Lysis Buf.2 (4X) 2 mL	12RA

Results :

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

[Apo C3 ] ng/ml	Conc. (ng/ml)	DF %	DF Accepted Range
Std 1	1,1	82	80 -119
Std 2	2,43	166	147-221
Std 3	5,34	344	295-443
Std 4	11,74	746	613-919
Std 5	25,83	1565	1327-1990
Std 6	56,82	3183	2740-4110
Std 7	125	3409	3183-4775



HTRF (h/m) HTT Mutant Kit - 500 pts - Lot n° : 03C

Accepted

Rejected

Quality control date : 2024-03-25	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.