

HTRF® Research Reagents

HTRF (h/m) HTT Mutant Kit - 10K pts

Product Number : 64HTTMPEH Lot Number : 02A

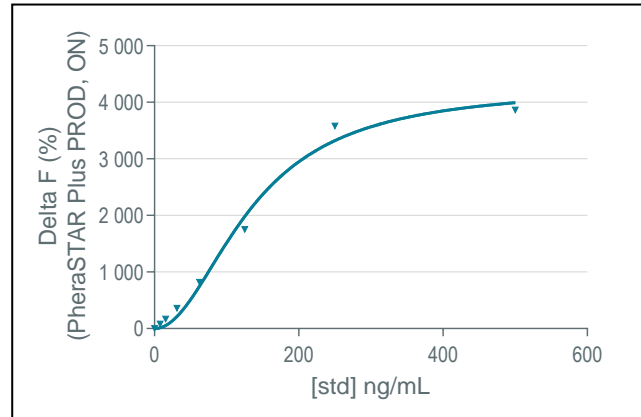
Expiration date : 2025-11-28 Document version : 1

Reagents :		Lot n°
HTRF Detect. Buf. 15 - 50 mL		06A
HTRF (h/m) HTT Mutant Kit - 10K pts d2 antibody		02A
HTRF (h/m) HTT Mutant Kit - 10K pts Tb Cryptate antibody		02A
HTRF (h/m) HTT Mutant Kit - modop		1
HTRF P-T prot. - Lysis Buf.2 (4X) 130 mL		12A

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	83%	80 -119
Std 2	2,43	169%	147-221
Std 3	5,34	362%	295-443
Std 4	11,74	819%	613-919
Std 5	25,83	1754%	1327-1990
Std 6	56,82	3581%	2740-4110
Std 7	125	3865%	3183-4775



HTRF (h/m) HTT Mutant Kit - 10K pts - Lot n° : 02A

 Accepted

 Rejected

Quality control date : 2022-02-12	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.