

HTRF® Research Reagents

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

Product Number : 64HTTTPEH Lot Number : 04A

Expiration date : 2025-10-03 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 15 - 50 mL	06A
HTRF (h/m) HTT Total Kit - 10K pts d2 antibody	04A
HTRF (h/m) HTT Total Kit - 10K pts Tb Cryptate antibody	04A
HTRF (h/m) HTT Total Kit - Ctrl Lysate	02A
HTRF (h/m) HTT Total 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 : pas de courbe

	Found value	Specifications Rev 01
Negative	962	
Control Lysate	7708	
S/B	8	>7

HTRF (h/m) HTT Total Kit - 10K pts - Lot n° : 04A

 Accepted

 Rejected

Quality control date : 2022_11_23	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.