

HTRF® Research Reagents

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

Product Number : 64HTTTPEH Lot Number : 04B

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

Reagents :	Lot n°
HTRF Detect. Buf. 15 - 2 mL	09RA
HTRF (h/m) HTT Total Kit - 500 pts d2 antibody	04AR
HTRF (h/m) HTT Total Kit - 500 pts Tb Cryptate antibody	04AR
HTRF (h/m) HTT Total Kit - Ctrl Lysate	02RA
HTRF (h/m) HTT Total Kit - modop	1
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 : pas de courbe

Assay characteristics :

	Found value	Specifications Rev 01
Negative	962	742-1187
Control Lysate	7708	5390-9808
S/B	8	>7

HTRF (h/m) HTT Total kit - 500 pts - Lot n° : 03B

 Accepted

 Rejected

Quality control date : 2023-09-20	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.