

## HTRF® Research Reagents

**HTRF Tau aggreg. kit - 10K pts**

Product Number : 6FTAUPEH Lot Number : 12D

Expiration date : 2025-12-21 Document version : 1

Reagents :	Lot n°
HTRF Diluent 3 - 20 mL	13AR
HTRF Tau aggreg. kit - 10K pts d2 antibody	12RA
HTRF Tau aggreg. kit - 10 K pts Tb Cryptate antibody	12RA
HTRF Tau aggreg. Kit - Positive Control	06RA
HTRF P-T prot. - Lysis Buf.1 (4X) 130mL	34RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B

**Results :**

 Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model :

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	939	/
Control Lysate	7246	/
S/B	7,7	>5
	/	/

HTRF Tau aggreg. kit - 10 K pts - Lot n° : 12D

 Accepted

 Rejected

Quality control date : 2024-11-08	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.