

HTRF® Research Reagents

HTRF Tau aggreg. kit - 500 pts

Product Number : 6FTAUPEG Lot Number : 12L

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

Reagents :	Lot n°
HTRF Diluent 3 - 20 mL	13AR
HTRF Tau aggreg. kit - 500 pts d2 antibody	12RA
HTRF Tau aggreg. kit - 500 pts Tb Cryptate antibody	12RA
HTRF Tau aggreg. Kit - Positive Control	06RA
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B

Results :

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

Assay characteristics :

	Found value	Specifications Rev 01
Negative	563	/
Control Lysate	4575	/
S/B	8,1	>5
	/	/

HTRF Tau aggreg. kit - 500 pts - Lot n° : 12L

 Accepted

 Rejected

Quality control date : 2024-11-04	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.