



## Quality Control report

HTRF Tau aggreg. kit - 10K pts

( 6FTAUPEH )

Lot n°12B

Expiration date : 2025-05- 19

Reagents :	Lot n°
HTRF Diluent 3 - 20 ml	11A
HTRF Tau aggreg. kit - 10K pts modop	8
HTRF Tau aggreg. kit - 10K pts d2 antibody	12RA
HTRF Tau aggreg. kit - 10K pts Tb Cryptate antibody	12RA
HTRF Tau aggreg. kit - Positive Control	06RA
HTRF P-T prot.- lysis buf.1 (4X) 130 mL	34RA
HTRF P-T prot. Block. Reag.(100X) 2 ml	28A

**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 - 10K pts Lot n°12B

Accepted

Rejected