

HTRF® Research Reagents

HTRF Histone H3 total kit - 500 pts

Product Number : 64NH3PEG Lot Number : 08A

Expiration date : 14/09/2024 Document version : 1

Reagents :	Lot n°
HTRF Histone H3 total kit - 500 pts d2 antibody	08A
HTRF Histone H3 total kit - 500 pts Eu Cryptate antibody	05A
HTRF Histone H3 total kit - Ctrl lysate	10A
HTRF Histone H3 total kit- modop	4
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34A
HTRF P-T prot. - Block. reag.(100X) 2 mL	26A
HTRF P-T prot. - Detect. Buf. 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	470	/
Control Lysate	3393	/
S/B	7,2	>5
	/	/

HTRF Histone H3 total kit - 500 pts - Lot n° : 08A
 Accepted

 Rejected

Quality control date : 2022-05-10 Supervisor's initials : AC