

HTRF® Research Reagents

## HTRF Histone H3 total kit - 500 pts

Product Number : 64NH3PEG Lot Number : 08D

Expiration date : Document version : 1

Reagents :	Lot n°
HTRF Histone H3 total kit - 500 pts d2 antibody	08BR
HTRF Histone H3 total kit - 500 pts Eu Cryptate antibody	06CR
HTRF Histone H3 total kit - Ctrl lysate	11AR
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Detect. Buf. 2 mL	32B

**Results :**

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

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	429	/
Control Lysate	4589	/
S/B	10.7	>5
	/	/

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

 Accepted

 Rejected

Quality control date : 2023-09-14 Supervisor's initials : AC

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.