

HTRF® Research Reagents

HTRF H3K36Me2 cell. assay - 500 pts

Product Number : 62KD2PAE Lot Number : 05Q

Expiration date : 15/07/2024 Document version : 1

Reagents :	Lot n°
HTRF Cellular Histone - Detect. Buf. 6 mL	31A
HTRF H3K36Me2 cell. assay - 500 pts d2 antibody	05RD
HTRF H3K36Me2 cell. assay - 500 pts Eu Cryptate antibody	04RD
HTRF H3K36Me2 cell. assay - Ctrl lysate	09RA
HTRF Epig. Lysis Buf. B - (4X)- 500 pts	06RB
HTRF Epig. Lysis Buf. C - 500 pts	11A

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	467	/
Control Lysate	4145	/
S/B	8,8	>5
	/	/

HTRF H3K36Me2 cell. assay - 500 pts - Lot n° : 05Q

 Accepted

 Rejected

Quality control date : 2023-03-17	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.