

HTRF® Research Reagents

HTRF (h) Cyclin E1 Total Kit - 10K pts

Product Number : 64CYCE1TPEH Lot Number : 01A

Expiration date : 2026-03-13 Document version : 1

Reagents :	Lot n°
HTRF (h) Cyclin E1 Total Kit - 10K pts d2 antibody	01A
HTRF (h) Cyclin E1 Total Kit - 10K pts Eu Cryptate antibody	01A
HTRF (h) Cyclin E1 Tot Kit - Ctrl Lys.	01A
HTRF (h) Cyclin E1 Total Kit - modop	1
HTRF P-T prot. - Lysis Buf.2 (4X) 130 mL	12RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28A
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

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	466	/
Control Lysate	3737	/
S/B	8	>5
/	/	/

HTRF (h) Cyclin E1 Total Kit - 10K pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2023-04-04	Supervisor's initials : PR
-----------------------------------	----------------------------

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.