

HTRF® Research Reagents

HTRF (h/m) CDK1 p-Y15 Kit - 500 pts

Product Number : 64CDK1Y5PEG Lot Number : 01E

Expiration date : 2025-11-03 Document version : 1

Reagents :	Lot n°
HTRF (h/m) CDK1 p-Y15 Kit - 500 pts d2 antibody	01RA
HTRF (h/m) CDK1 p-Y15 Kit - 500 pts Eu Cryptate antibody	01RA
HTRF (h/m) CDK1 p-Y15 Kit - Ctrl Lysate	01RA
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28RA
HTRF P-T prot. - Detect. Buf. 2 mL	32B

Results :

 Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	512	/
Control Lysate	5882	/
S/B	11.5	>5
	/	/

HTRF (h/m) CDK1 p-Y15 Kit - 500 pts - Lot n° : 01E

 Accepted

 Rejected

Quality control date : 2023-01-12 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.