

## HTRF® Research Reagents

## HTRF Human AR-V7 Kit - 500 pts

Product Number : 64ARV7PEG

Lot Number : 02A

Expiration date : 2026-04-03

Document version : 1

Reagents :		Lot n°
HTRF Human AR-V7 Kit - 500 pts d2 antibody		02A
HTRF Human AR-V7 Kit - 500 pts Eu Cryptate antibody		02A
HTRF Human AR-V7 Kit - Ctrl Lysate		02A
HTRF P-T prot. - Lysis Buf.3 (4X) 2 mL		14B
HTRF P-T prot. - Block. reag.(100X) 0.3 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 : EnVision

Fitting model : /

## Assay characteristics :

	Found value	Specifications Rev 01
Negative	386	/
Control Lysate	2970	/
S/B	7.7	>5

HTRF Human AR-V7 Kit - 500 pts - Lot n° : 02A

☒ Accepted☐ Rejected

Quality control date : 2024-10-04

Supervisor's initials : ml

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.