

**HTRF® Research Reagents**
**HTRF Human AR-V7 Kit - 10K pts**

Product Number : 64ARV7PEH Lot Number : 03A

Expiration date : 2026-04-03 Document version : 1

Reagents :	Lot n°
HTRF Human AR-V7 Kit - 10K pts d2 antibody	03A
HTRF Human AR-V7 Kit - 10K pts Eu Cryptate antibody	03A
HTRF Human AR-V7 Kit - Ctrl Lysate	02A
HTRF P-T prot. - Lysis Buf.3 (4X) 130 mL	14RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
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	494	/
Control Lysate	2934	/
S/B	5.9	>5
	/	/

HTRF Human AR-V7 Kit - 10K pts - Lot n° : 03A

☒ Accepted

☐ Rejected

Quality control date : 2024-12-03 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.