

## HTRF® Research Reagents

# HTRF Andro Receptor Kit - 10K pts

Product Number : 64ANDRPEH Lot Number : 05A

Expiration date : 2026-07-05 Document version : 1

Reagents :	Lot n°
HTRF Andro Receptor Kit - 10K pts d2 antibody	05A
HTRF Andro Receptor Kit - 10K pts Eu Cryptate antibody	05A
HTRF Andro Receptor Kit - Ctrl lysate	04AR
HTRF P-T prot. - Lysis Buf.3 (4X) 130 mL	14RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	29A
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

**Results :**

 Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : No curve

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	1182	/
Control Lysate	11180	/
S/B	11,1	>5

HTRF Andro Receptor Kit - 10K pts - Lot n° : 05A

 Accepted

 Rejected

Quality control date : 2024-12-27	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.