

## HTRF® Research Reagents

# HTRF GAPDH Housekeeping kit - 500 pts

Product Number : 64GAPDHPEG Lot Number : 07D

Expiration date : 2027-11-07 Document version : 1

Reagents :	Lot n°
HTRF Diluent 11 - 10 mL	03C
HTRF GAPDH Housekeeping kit - 500 pts d2 antibody	07A
HTRF GAPDH Housekeeping kit - 500 pts Eu Cryptate antibody	07A
HTRF GAPDH House. kit - Ctrl lysate	07A
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	29A
HTRF P-T prot. - Detect. Buf. 2 mL	33A

**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	509	/
Control Lysate	2763	/
S/B	5,4	>5

HTRF GAPDH Housekeeping kit - 500 pts - Lot n° : 07D

 Accepted

 Rejected

Quality control date : 2025-03-07	Supervisor's initials : LM
-----------------------------------	----------------------------

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.