

HTRF® Research Reagents

## HTRF Active GLP1 kit - 500 pts

Product Number : 62GLPPEG Lot Number : 11E

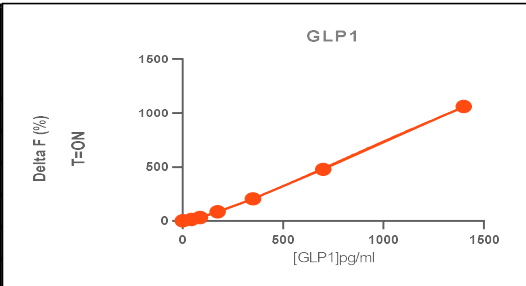
Expiration date : 2027-07-04 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 5 - 7 mL	07B
HTRF Active GLP1 kit - Std	07A
HTRF Diluent 4 - 10 mL	10A
HTRF Active GLP1 kit - 500 pts d2 antibody	11B
HTRF Active GLP1 kit - 500 pts Tb Cryptate antibody	11B

## Results :

Incubation : Overnight at RT  
Plate : d  
Reader : Pherastar Plus  
Fitting model : 4 Parameter Logistic (4PL)

GLP1	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	43,75	13	6-13
Std 2	87,5	32	14-37
Std 3	175	86	42-93
Std 4	350	205	125-239
Std 5	700	477	335-619
Std 6	1400	1061	833-1456



## Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	13 pg/ml	<=25 pg/ml

HTRF Active GLP1 kit - 500 pts - Lot n° : 11E

☒ Accepted☐ Rejected

Quality control date : 2024-11-05	Supervisor's initials : MP
-----------------------------------	----------------------------

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.