

HTRF® Research Reagents

HTRF Active GLP1 kit - 500 pts

Product Number : 62HGRZBPEG Lot Number : 11B

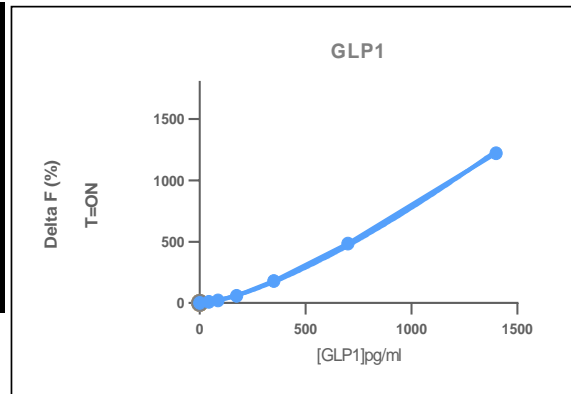
Expiration date : 24/07/2024 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 5 - 7 mL	07RA
HTRF Active GLP1 kit Std	06RA
HTRF Diluent 4 - 10 ml	08RA
HTRF Active GLP1 kit - 500 pts d2 antibody	11RA
HTRF Active GLP1 kit - 500 pts Tb Cryptate antibody	11RA
HTRF Active GLP1 kit - modop	4

Results :

Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µl  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL)

GLP1	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	43,75	9%	6-13
Std 2	87,5	22%	14-37
Std 3	175	59%	42-93
Std 4	350	179%	125-239
Std 5	700	485%	335-619
Std 6	1400	1220%	833-1456



Assay characteristics :

	Found value	Specifications
Detection limite	20,7	<=25 pg/ml

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

Accepted

Rejected

Quality control date : 2023-09-01	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