

HTRF® Research Reagents

HTRF Active GLP1 kit - 500 pts

Product Number : 62GLPPEG Lot Number : 10A

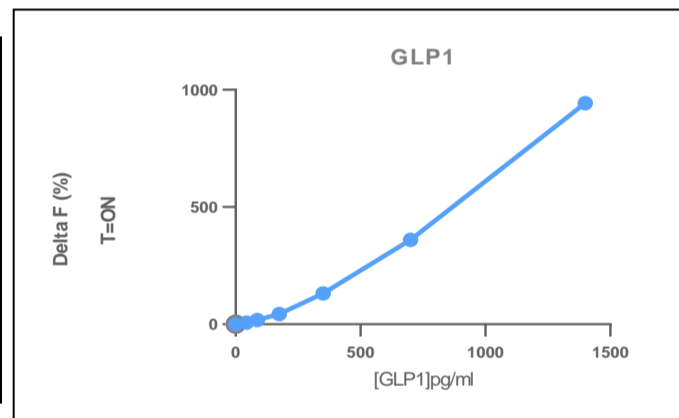
Expiration date : 2026-07-24 Document version : 1

Reagents :		Lot n°
HTRF Detect. Buf. 5 - 7 mL		08AR
HTRF Active GLP1 kit - Std		06RA
HTRF Diluent 4 - 10 mL		09AR
HTRF Active GLP1 kit - 500 pts d2 antibody		10RB
HTRF Active GLP1 kit - 500 pts Tb Cryptate antibody		10RB

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	10%	6-13
Std 2	87,5	32%	14-37
Std 3	175	64%	42-93
Std 4	350	187%	125-239
Std 5	700	357%	335-619
Std 6	1400	1027%	833-1456



Assay characteristics :

	Found value	Specifications
Detection limite	17,7	<=25 pg/ml

HTRF Active GLP1 kit - 500 pts - Lot n° : 10A

Accepted

Rejected

Quality control date : 2024-02-02	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.