

HTRF® Research Reagents

HTRF Active GLP1 kit - 500 pts

Product Number : 62GLPPEG Lot Number : 11C

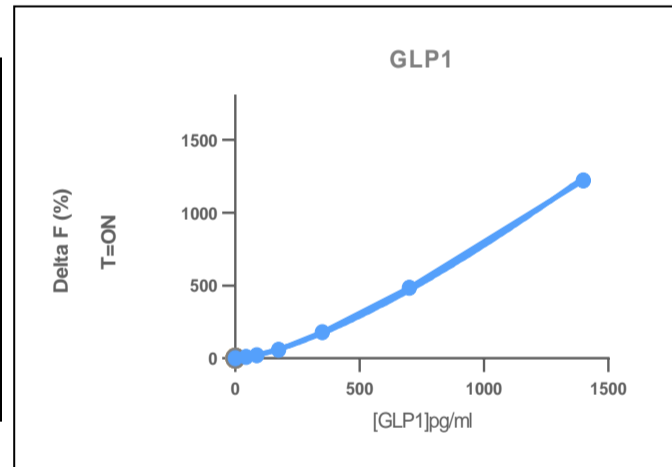
Expiration date : 2025-03-31 Document version : 1

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

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 Rev 01
Detection limite	20,7	<=25 pg/ml

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

Accepted

Rejected

Quality control date : 2024-02-05	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.