

HTRF® Research Reagents

HTRF Active GLP1 kit - 10K pts

Product Number : 62GLPPEH Lot Number : 10D

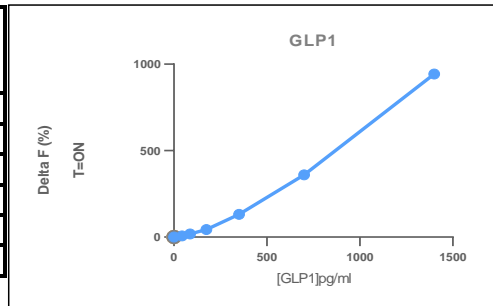
Expiration date : 2024-07-24 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 5 - 105 mL	07A
HTRF Active GLP1 kit - 10K pts d2 antibody	10RA
HTRF Active GLP1 kit - 10K pts Tb Cryptate antibody	10RA
HTRF Active GLP1 kit - Std	06RA
HTRF Diluent 4 - 10 mL	08RA
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	7%	6-13
Std 2	87,5	19%	14-37
Std 3	175	44%	42-93
Std 4	350	132%	125-239
Std 5	700	361%	335-619
Std 6	1400	943%	833-1456


Assay characteristics :

	Found value	Specifications
Detection limite	23,6	<=25 pg/ml

HTRF Active GLP1 kit - 10K pts - Lot n° : 10D

 Accepted

 Rejected

Quality control date : 2023-11-09	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.