

HTRF® Research Reagents

HTRF Glucagon kit - 500 pts

Product Number : 62CGLPEG Lot Number : 06H

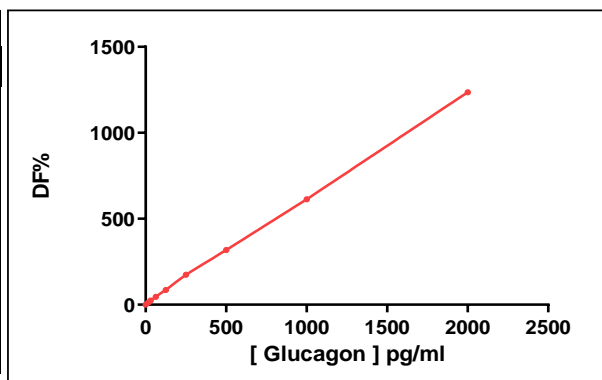
Expiration date : 23/12/2024 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 6 - 7 mL	10RB
HTRF Glucagon kit - 500 pts d2 antibody	06RA
HTRF Glucagon kit - 500 pts Tb Cryptate antibody	06RA
HTRF Glucagon kit - modop	4
HTRF Glucagon kit - Std	11C

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : Linear regression

Glucagon] (pg/ml)	DF %	Diluent #5	
		Range	
15,6	12%	8	13
31,3	24%	18	25
62,5	45%	36	49
125	86%	68	92
250	174%	145	183
500	318%	297	355
1000	613%	595	701
2000	1236%	1234	1531



HTRF Glucagon kit - 500 pts - Lot n° : 06H

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 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.