

HTRF® Research Reagents

HTRF Glucagon kit - 500 pts

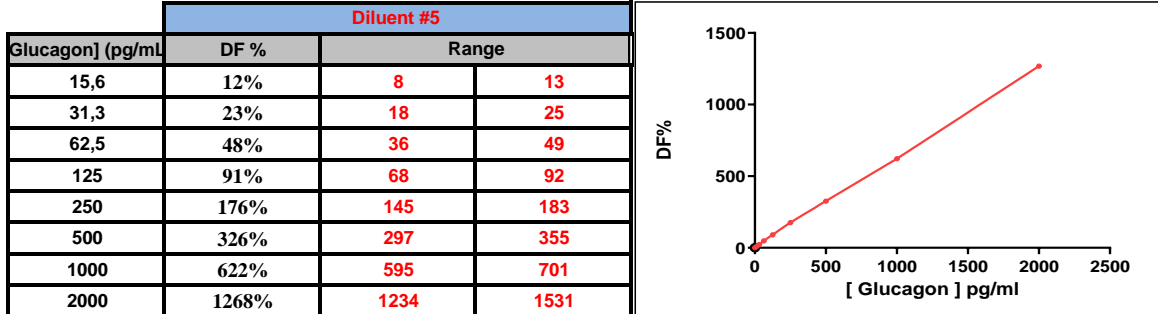
Product Number : 62CGLPEG Lot Number : 01B

Expiration date : 2026-12-20 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		01C
HTRF Glucagon kit - 500 pts Tb Cryptate antibody		01C
HTRF Glucagon kit - Std		12AR

Results :

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



Assay characteristics :

	Found value	Specifications
Detection limite	5,3 pg/ml	< 7 pg/ml

HTRF Glucagon kit - 500 pts - Lot n° : 01B

Accepted

Rejected

Quality control date : 2025-02-28	Supervisor's initials : PL
-----------------------------------	----------------------------

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.