

HTRF® Research Reagents

### HTRF Insulin High Range kit - 500 pts

Product Number : 62IN1PEG Lot Number : 09C

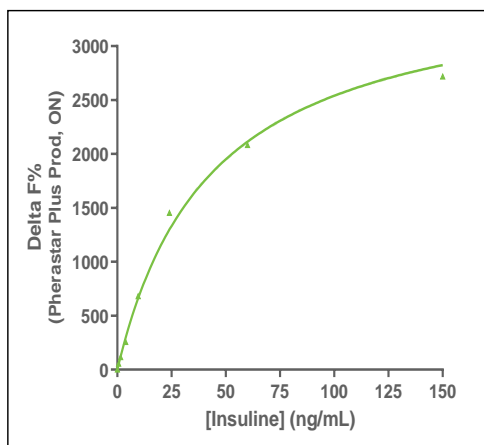
Expiration date : 2026-06-02 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 7 - 55 mL	13RA
HTRF Insulin High Range kit - 500 pts Eu Cryptate antibody	09RA
HTRF Insulin High Range kit - 500 pts XL665 antibody	09RA
HTRF Insulin High Range kit - modop	5
HTRF Insulin kit - Std	16RA

**Results :**

Incubation : Overnight at RT  
 Plate : 384 wells regular white plate 80µL  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL)

Insulin standards	[Insulin] (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0	0-0
Std 1	0,6144	56%	35-77
Std 2	1,536	116%	80-197
Std 3	3,84	256%	200-452
Std 4	906	682%	500-930
Std 5	24	1454%	1000-1840
Std 6	60	2083%	2000-2650
Std 7	150	2718%	2700-3513



**Assay characteristics :**

	Found value	Specifications Rev 01
Detection limit	0.044 ng/ml	0.02-0.06 ng/ml

HTRF Insulin High Range kit - 500 pts - Lot n° : 09C

Accepted

Rejected

Quality control date : 2023-11-13	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.