

HTRF® Research Reagents

HTRF Insulin High Range kit - 500 tests

Product Number : 62IN1PEG Lot Number : 09B

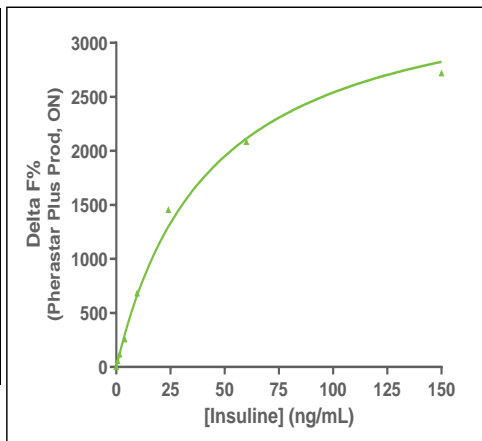
Expiration date : 02/06/2024 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21A
HTRF Detect. Buf. 3 - 1.5 mL		13RA
HTRF Human Granzyme B kit - Std		16A
HTRF Human Granzyme B kit - 500 pts d2 antibody		09A
HTRF Human Granzyme B kit - 500 pts Eu Cryptate antibody		09A
HTRF Human Granzyme B kit - modop		5

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 - £ 09B

 Accepted

 Rejected

Quality control date : 2023-09-11	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