

HTRF® Research Reagents

HTRF Insulin High Range kit - 10K pts

Product Number : 62IN1PEH Lot Number : 08B

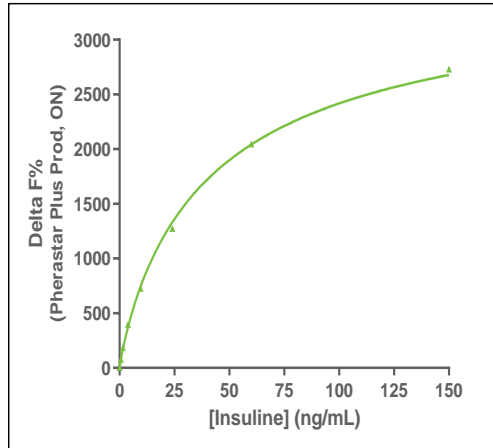
Expiration date : 2026-01-25 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 100 mL	21RA
HTRF Detect. Buf. 7 - 130 mL	13RA
HTRF Insulin High Range kit - 10K pts Eu Cryptate antibody	08RA
HTRF Insulin High Range kit - 10K pts XL665 antibody	08RA
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	76%	35-77
Std 2	1,536	183%	80-197
Std 3	3,84	394%	200-452
Std 4	906	724%	500-930
Std 5	24	1271%	1000-1840
Std 6	60	2043%	2000-2650
Std 7	150	2729%	2700-3513



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	0.032 ng/ml	0.02-0.06 ng/ml

HTRF Insulin High Range kit - 10K pts - Lot n° : 08B

Accepted

Rejected

Quality control date : 2023-11-30	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.