

HTRF® Research Reagents

HTRF Insulin High Range kit - 500 pts

Product Number : 62IN1PEG Lot Number : 16A

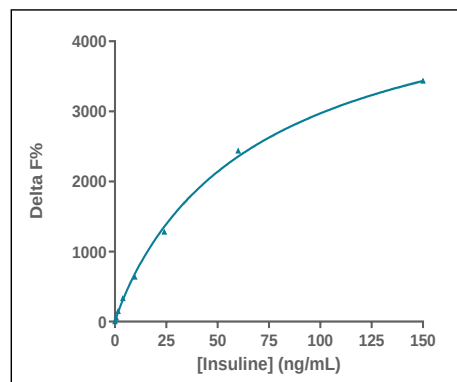
Expiration date : 2028-05-26 Document version :

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	22A
HTRF Detect. Buf. 7 - 55 mL	14AR
HTRF Insulin High Range kit - 500 pts Eu Cryptate antibody	16A
HTRF Insulin High Range kit - 500 pts XL665 antibody	16A
HTRF Insulin kit - Std	18A

Results :

Incubation : Overnight at RT
Plate : 384 wells regular white plate 80µL
Reader : EnVision
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	54%	35-77
Std 2	1,536	146%	80-197
Std 3	3,84	331%	200-452
Std 4	906	637%	500-930
Std 5	24	1280%	1000-1840
Std 6	60	2436%	2000-2650
Std 7	150	3434%	2700-3513



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	0.04 ng/ml	0.02-0.06 ng/ml

HTRF Insulin High Range kit - 500 pts - Lot n° : 16A

☒ Accepted☐ Rejected

Quality control date : 2025-07-02	Supervisor's initials : FL
-----------------------------------	----------------------------

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.