

HTRF® Research Reagents

HTRF Insulin Mouse Serum kit - 5x200 pts

Product Number : 62IN3PEB Lot Number : 03M

Expiration date : 2024-05-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 6 - 6 mL	15RA
HTRF Detect. Buf. 6 - 3 mL	12RA
HTRF Insulin Mouse Serum kit - 200 pts modop	5
HTRF Insulin Mouse Serum kit - 200 pts Tb Cryptate antibody	03RA
HTRF Insulin Mouse Serum kit - 200 pts XL665 antibody	03RA
HTRF Insulin kit - Std	16RA

Results :

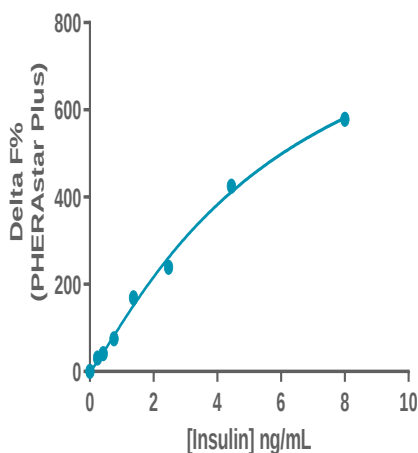
Incubation : Overnight at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar Plus

Fitting model : 4 Parameter Logistic (4PL)

Insulin standards	[Insulin] (ng/ml)	Delta F %	Delta F % Accepted Range Rev 2
Std 0	0	0	0-0
Std 1	0,24	31	/
Std 2	0,42	41	33 - 47
Std 3	0,76	75	58 - 96
Std 4	1,37	169	/
Std 5	2,47	239	180 - 341
Std 6	4,44	425	/
Std 7	8	578	545 - 758



Assay characteristics :

	Found value	Specifications
Detection limite	0,031	0.055 ng/ml

HTRF Insulin Mouse Serum kit - 5x200 pts - Lot n° : 03M

☒ Accepted☐ Rejected

Quality control date :	2023-11-10	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.