

HTRF® Research Reagents

HTRF Insulin Mouse Serum kit - 5x200 pts

Product Number : 62IN3PEB

Lot Number : 03P

Expiration date : 2026-05-16

Document version : 1

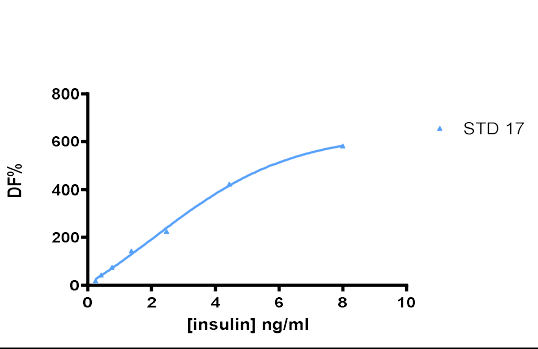
Reagents :	Lot n°
HTRF Diluent 6 - 6 mL	16AR
HTRF Detect. Buf. 6 - 3 mL	13AR
HTRF Insulin Mouse Serum kit - 200 pts Tb Cryptate antibody	03RA
HTRF Insulin Mouse Serum kit - 200 pts XL665 antibody	03RA
HTRF Insulin kit - Std	17A

Results :

Incubation : Overnight at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : 4 Parameter Logistic (4PL)

Assay characteristics :

Insulin standards	[Insulin] (ng/ml)	Delta F %	Delta F % Accepted Range Rev 2
Std 0	0	0	0-0
Std 1	0.24	18	13 - 25
Std 2	0.42	42	33 - 47
Std 3	0.76	74	58 - 96
Std 4	1.37	142	96 - 180
Std 5	2.47	225	180 - 341
Std 6	4.44	422	341 - 545
Std 7	8	582	545 - 758



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

☒ Accepted☐ Rejected

Quality control date : 2024-01-12

Supervisor's initials : AC

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.