

HTRF® Research Reagents

# HTRF Insulin High Range kit - 10K pts

Product Number : 62IN1PEH Lot Number : 14A

Expiration date : 2027-01-03 Document version : 1

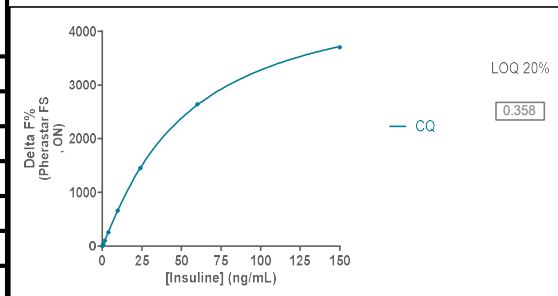
| Reagents :   | Lot n° |
|--|--------|
| HTRF Diluent 5 (5X) - 100 mL                               | 21RA   |
| HTRF Detect. Buf. 7 - 130 mL                               | 13RA   |
| HTRF Insulin High Range kit - 10K pts Eu Cryptate antibody | 14A    |
| HTRF Insulin High Range kit - 10K pts XL665 antibody       | 14A    |
| HTRF Insulin kit - Std                                     | 17A    |

**Results :**

Incubation : Overnight at RT  
 Plate : 384 wells regular white plate 80µL  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL)

**Assay characteristics :**

| Insulin standards | [Insulin] (ng/ml) | Delta F % | Delta F% Accepted Range |
|-------------------|-------------------|-----------|-------------------------|
| Std 0             | 0                 | 0         | 0-0                     |
| Std 1             | 0.6144            | 36%       | 35-77                   |
| Std 2             | 1.536             | 100%      | 80-197                  |
| Std 3             | 3.84              | 255%      | 200-452                 |
| Std 4             | 906               | 658%      | 500-930                 |
| Std 5             | 24                | 1452%     | 1000-1840               |
| Std 6             | 60                | 2640%     | 2000-2650               |
| Std 7             | 150               | 3703%     | 2700-3513               |



HTRF Insulin High Range kit - 10K pts - Lot n° : 14A

Accepted

Rejected

Quality control date : 2024-07-23

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.