

HTRF® Research Reagents

HTRF HIV p24 kit - 10K pts

Product Number : 64P24PEH Lot Number : 11A

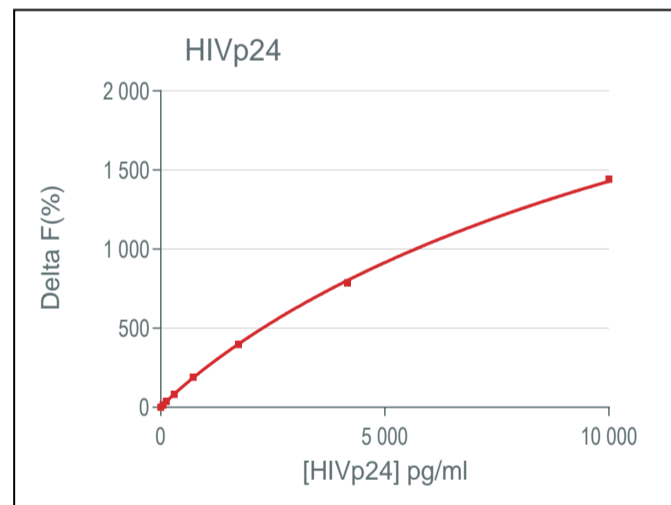
Expiration date : 2026-06-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF HIV p24 kit - 10K pts d2 antibody	11A
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	11A
HTRF HIV p24 kit - Std	05A

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL)

HIV p24	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 1	52	16%	oct-23
Std 2	126	38%	22 -55
Std 3	301	83%	56 - 124
Std 4	723	190%	135 - 273
Std 5	1736	398%	303 - 553
Std 6	4167	788%	644 - 1003
Std 7	10000	1442%	1023 - 1524



HTRF HIV p24 kit - 10K pts - Lot n° : 11A

Accepted

Rejected

Quality control date : 2024-07-12	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.