

HTRF® Research Reagents

# HTRF HIV p24 kit - 500 pts

Product Number : 64P24PEG Lot Number : 02F

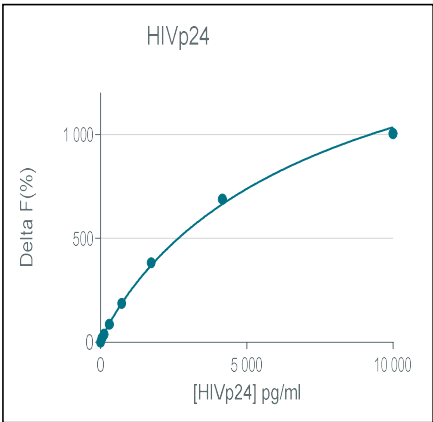
Expiration date : 2025-05-09 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF HIV p24 kit - Std	04B
HTRF HIV p24 kit - 500 pts d2 antibody	02RA
HTRF HIV p24 kit - 500 pts Eu Cryptate antibody	02RA

**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	20%	14-21
Std 2	126	38%	28-41
Std 3	301	86%	61-92
Std 4	723	187%	145-217
Std 5	1736	383%	307-460
Std 6	4167	689%	600-900
Std 7	10000	1005%	971-1456



HTRF HIV p24 kit - 500 pts - Lot n° : 02F  Accepted  Rejected

Quality control date : 2023-12-21 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.