

HTRF® Research Reagents

HTRF HIV p24 kit - 500 pts

Product Number : 64P24PEG Lot Number : 02D

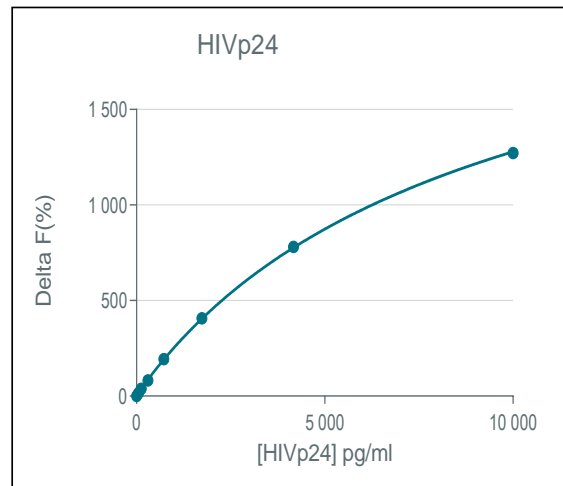
Expiration date : 2025-05-09 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21RA
HTRF Detect. Buf. 3 - 1.5 mL		28RA
HTRF HIV p24 kit - Std		03RA
HTRF HIV p24 kit - 500 pts d2 antibody		02RA
HTRF HIV p24 kit - 500 pts Eu Cryptate antibody		02RA
HTRF HIV p24 kit - modop		2

Results :

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

HMGB1	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 1	52	17%	14-21
Std 2	126	38%	28-41
Std 3	301	86%	61-92
Std 4	723	201%	145-217
Std 5	1736	422%	307-460
Std 6	4167	814%	600-900
Std 7	10000	1323%	971-1456



HTRF HIV p24 kit - 500 pts - Lot n° : 02D

Accepted

Rejected

Quality control date : 2023-10-18	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.