

HTRF® Research Reagents

HTRF AAV5 Capsid Kit - 500 pts

Product Number : 64AAV5PEG Lot Number : 01A

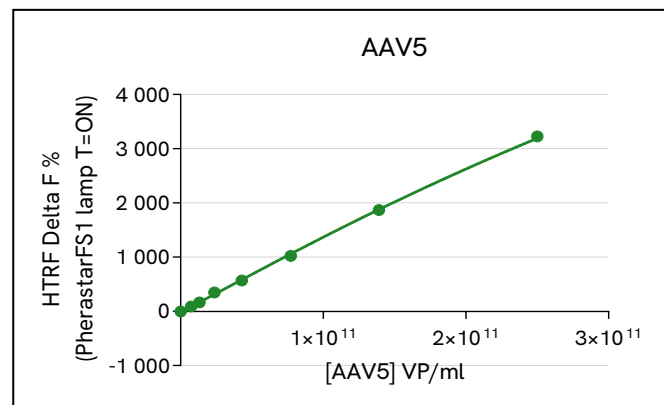
Expiration date : 2026-06-26 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF AAV5 Capsid Kit - Std	01A
HTRF AAV5 Capsid Kit - 500 pts d2 antibody	01A
HTRF AAV5 Capsid Kit - 500 pts Biotin antibody	01A
HTRF AAV5 Capsid Kit - 500 pts Eu Cryptate Streptavidin	01A
HTRF P-T prot. - Lysis Buf.3 (4X) 2 mL	14A

Results :

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

	Conc (VP/ml)	Delta F %	Delta F% Accepted Range
Std 0	0,00E+00	0%	0-0
Std 1	7,35E+09	89%	31-147
Std 2	1,32E+10	165%	67-263
Std 3	2,38E+10	346%	132-560
Std 4	4,29E+10	568%	372-764
Std 5	7,72E+10	1024%	650-1398
Std 6	1,39E+11	1868%	1246-2491
Std 7	2,50E+11	3226%	1862-4589



HTRF AAV5 Capsid Kit - 500 pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2023-12-22	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.