

HTRF® Research Reagents

HTRF AAV9 Capsid Kit - 500 pts

Product Number : 64AAV9PEG Lot Number : 01B

Expiration date : 2026-06-19 Document version : 1

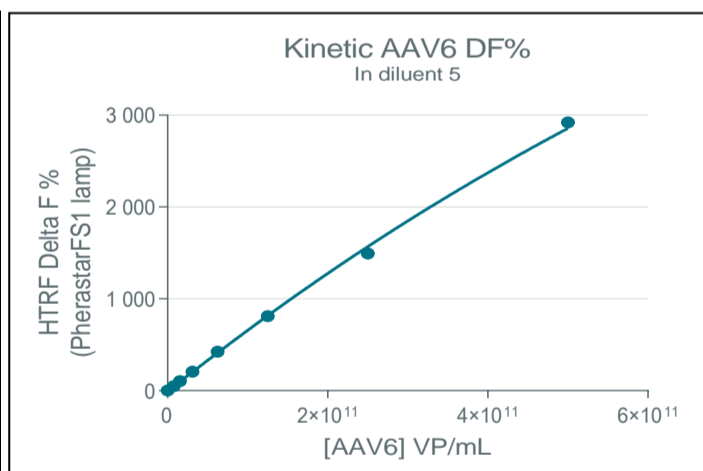
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF AAV9 Capsid Kit - Std	01A
HTRF AAV9 Capsid Kit - 500 pts d2 antibody	01A
HTRF AAV9 Capsid Kit - 500 pts Eu Cryptate	01A
HTRF P-T prot. - Lysis Buf.3 (4X) 2 mL	14B

Results :

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

Assay characteristics :

	Conc (VP/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	7812500000	48%	28-68
Std 2	15625000000	103%	67-138
Std 3	31250000000	206%	144-268
Std 4	62500000000	424%	297-551
Std 5	1,25E+11	810%	591-1029
Std 6	2,5E+11	1494%	1279-1709
Std 7	5E+11	2920%	2322-3517



HTRF AAV9 Capsid Kit - 500 pts - Lot n° : 01B

Accepted

Rejected

Quality control date : 2024-08-23	Supervisor's initials : PL
-----------------------------------	----------------------------

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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.