



AAVPrime™ AAV Serotype Testing Kit

Cat. No. AA320

AAVPrime™ AAV Serotype Testing Kit Plus

Cat. No. AA321

Datasheet

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, MD 20850 USA

Tel: 301-762-0888, Fax: 301-762-3888

Email: inquiry@genecopoeia.com

Website: www.genecopoeia.com

I. Introduction

GeneCopoeia's AAVPrime™ Adeno-associated virus (AAV) Serotype Testing Kit (Cat. No. AA320) contains 9 premade GFP-expressing AAV in serotypes 1, 2, 3, 6, 7, 8, 9, DJ and DJ/8. AAVPrime™ Adeno-associated virus (AAV) Serotype Testing Kit Plus (Cat. No. AA321) contains 12 premade GFP-expressing AAV in serotypes 1, 2, 3, 4, 5, 6, 7, 8, 9, 10, DJ and DJ/8. All AAV particles are offered at 5×10^{12} GC/ml, in 25 μ l stocks, except serotypes 4 and 10 are offered at 5×10^{11} - 10^{12} GC/ml. These kits are ideal for testing the infection efficiency of AAV serotypes in different cell types.

II. Contents and Storage

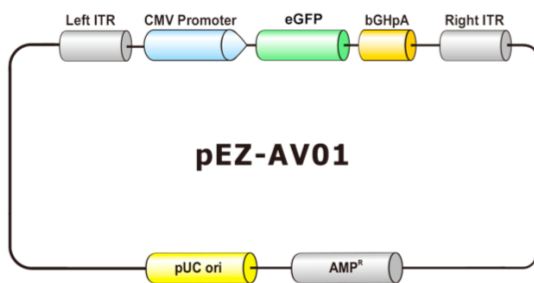
AAV stocks are supplied in liquid form at the indicated titer in PBS storage solution with 5% glycerol. Store at -80 °C. If desired, aliquot viral stock upon arrival, and store the aliquots at -80 °C immediately. Avoid repeated freeze and thaw.

Catalog	Products	Titer	Quantity
AA601	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 1. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AA001	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 2. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AA801	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 3. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AB001*	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 4. Promoter: CMV.	$\geq 5 \times 10^{11}$ GC/ml	25 μ l
AB201*	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 5. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AB401	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 6. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AB601	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 7. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AB801	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 8. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AC001	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 9. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l

AC201*	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: 10. Promoter: CMV.	$\geq 5 \times 10^{11}$ GC/ml	25 μ l
AA201	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: DJ. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l
AA401	eGFP AAVPrime™ Purified Adeno-associated Viral Particles. Serotype: DJ/8. Promoter: CMV.	$\geq 5 \times 10^{12}$ GC/ml	25 μ l

* AB001, AB201, and AC201 are not included in AAVPrime™ AAV Serotype Testing Kit (AA320).

III. Plasmid Map



IV. Quality Control

Every lot is tested for performance consistency, titer, sterility and purity.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at (301)-762-0888.