

Enhanced Safety • High Titers

AAVPrime™ Adeno-associated Viral (AAV) Particles

For *in vitro* and *in vivo* studies



Advantages

- Titer of purified AAV particles can be up to 10^{14} GC/ml
- All serotypes available - allows tissue selectivity
- Purified particles for *in vivo* animal studies
- Low toxicity and minimal host immune response
- Custom packaging service for ORF, shRNA, CRISPR and more

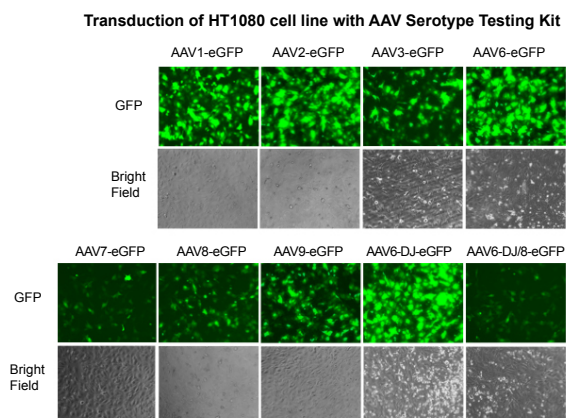
Primary Target Tissues

Serotype	Retina	Neurons	Brain	Lung	Heart	Liver	Muscle	Kidney	Pancreas
AAV-1		√			√		√		√
AAV-2	√	√	√			√	√	√	
AAV-3	√			√		√	√		
AAV-4	√	√	√				√		
AAV-5	√	√		√					
AAV-6				√	√	√	√		
AAV-7	√	√				√	√		√
AAV-8	√		√			√	√		
AAV-9			√	√	√	√	√	√	√
AAV-10		√		√	√	√	√		
AAV-DJ	Efficiently transduces a wide variety of cell types <i>in vitro</i>								
AAV-DJ/8	A variant of AAV-DJ that permits infection of liver as well as other tissues <i>in vivo</i>								

AAV Serotype Testing Kit

GeneCopoeia's AAVPrime™ Adeno-associated virus (AAV) Serotype Testing Kit contains 9 premade GFP-expressing AAV in serotypes 1, 2, 3, 6, 7, 8, 9, DJ and DJ/8. The kit can be used to determine the ideal AAV serotype for the infection of different tissue or cell types.

Kit	AAV
Components	vectors of 9 common serotypes
Size	9 x 25 µl
Titer	5×10^{12} GC/ml



AAV of different serotypes carrying GFP were used to transduce HT1080 cell line at MOI 20,000 as shown in the figure above. Fluorescent images (GFP) show that cells are transduced with AAV-GFP vectors from the AAV Serotype Testing Kit.

AAVPrime™ Adeno-associated Viral (AAV) Particles

Applications: For *in vivo* and *in vitro* studies of gene overexpression, knockdown and knockout. Type of custom AAV service available: ORF cDNA, shRNA, CRISPR sgRNA and SaCas9. Control AAV particles such as eGFP, mCherry, LacZ, luciferase, etc. in various serotypes are also available.

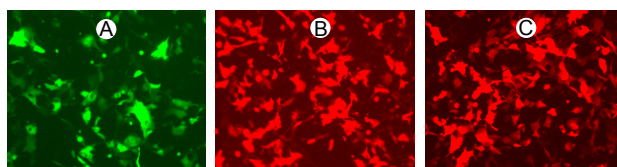


Figure 1. HT1080 cells in 24-well plates were transduced with 0.5 µL of standard AAV particles (MOI=200) expressing either eGFP (A), RFP (B), or mCherry (C). Cells were visualized with a fluorescence microscope 48 hours post-transduction (Exposure time: 400 ms).

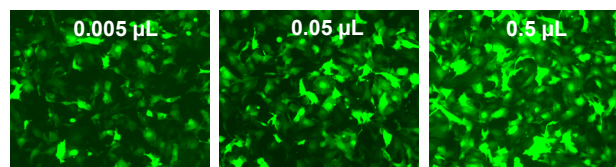


Figure 2. HT1080 cells in 24-well plates were transduced with the indicated amounts of purified AAV particles (MOI=100, 1000, 10000) expressing eGFP. Cells were visualized with a fluorescence microscope 72 hours post-transduction (Exposure time: 400 ms).

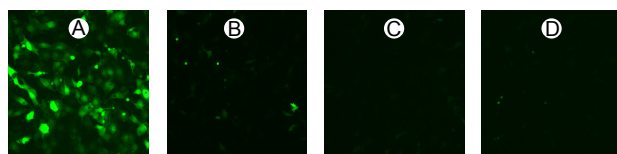


Figure 3. HT1080 cells in 24-well plates were co-transduced standard AAV particles AAV-2-eGFP expressing eGFP and AAV-2-shRNA knocking down eGFP with series of MOI ratios, 3000:0(A); 3000:1500(B); 3000:3000(C) and 3000:6000(D). Cells were visualized with a fluorescence microscope 48 hours post-transduction (Exposure time: 400 ms).

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA



Email
Tel
Toll free
Fax
Website

inquiry@genecopoeia.com
+1 (301) 762-0888
+1 (866) 360-9531
+1 (301) 762-3888
www.genecopoeia.com