

## AAVPrime™ AAV Serotype Testing Kit

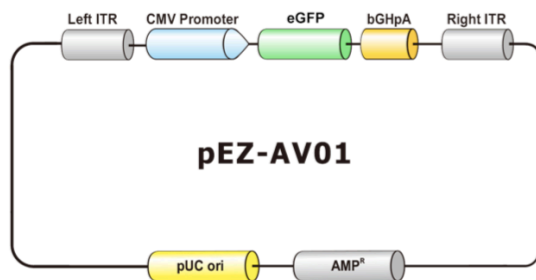
### I. Introduction

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. All AAV particles are offered in  $5 \times 10^{12}$  GC/ml, 25  $\mu$ l stocks.

### 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.

### 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.

## AAVPrime™ AAV Serotype Testing Kit Data Sheet

GeneCopoeia, Inc.

9620 Medical Center Drive, #101  
Rockville, Maryland 20850

Tel: 301-762-0888 Fax: 301-762-8333

Email: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

Web: [www.genecopoeia.com](http://www.genecopoeia.com)

GeneCopoeia Products are for Research Use Only

Copyright © 2018 GeneCopoeia, Inc.

Trademarks: GeneCopoeia™, AAVPrime™, (GeneCopoeia Inc.);

AAV-082118