

## CRE AAVPrime™ Adeno-Associated Virus - Serotype DJ (AAV-DJ) Particles

Cat. No. AA231-025, AA231-100

(Old Cat. No. AA231, AA232)

AAV-DJ infects a broad range of cell lines including both dividing and non-dividing mammalian cells and has been used successfully to express human and non-human proteins. AAV is not associated with any human disease, ensuring highly-efficient and safe results.

### Description

Gene: CRE  
Promoter: CMV  
Tag: CRE  
Reporter: N/A  
Additional note: N/A

### Contents and storage

Provided as 1 vial of 25 µl or 4 vials of 25 µl of purified AAV-DJ particles with titers of  $5 \times 10^{12}$  Genome Copies/ml (GC/ml).

AAV-DJ particles are shipped on dry ice and **must be stored at –80°C immediately upon receipt**. Avoid repeated freeze-thaw cycles as this will reduce titers. Divide into useful aliquots if necessary and store at –80°C.

### Viral titer

The physical copy number of the viral genome was determined by qPCR.

### Overview of production

Identity of AAV transfer vector: pEZ-AV01

The AAV particles were generated by following a standardized protocol using highly purified plasmids and EndoFectin-AAV™ reagents.

The AAV transfer vector was co-transfected into HEK293 cells with AAV-RC (serotype DJ) and Helper plasmids for AAV production.

### Quality control

The AAV expression construct was validated by full-length sequencing, restriction enzyme digestion and PCR-size validation using gene-specific and vector-specific primers. Product is confirmed free of bacteria, fungi and common *Mycoplasma* contamination.

GeneCopoeia, Inc.  
9620 Medical Center Drive, Suite 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)