

GLuc AAVPrime™ Adeno-Associated Virus - Serotype 2 (AAV-2) Particles

Cat. No. AA091-025, AA091-100

(Old Cat. No. AA091, AA092)

AAV-2 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: N/A
Promoter: CMV
Tag: N/A
Reporter: GLuc
Additional note: N/A

Contents and storage

Provided as 1 vial of 25 µl or 4 vials of 25 µl of purified AAV-2 particles with titers of 5×10^{12} Genome Copies/ml (GC/ml).

AAV-2 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 2) 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
Web: www.genecopoeia.com