

## ***iLite*<sup>®</sup> AAV8 Packaging Assay Ready Cells**

REF: BM6008

*For research use only. Not for use in diagnostic procedures.*

### **DESCRIPTION**

*iLite*<sup>®</sup> AAV8 Packaging Assay Ready Cells are based on the human embryonic kidney 293 (HEK293) cell line and have been genetically engineered and optimized to produce AAV vectors of serotype 8.

### **CONTENT**

>250 µL of Assay Ready Cells suspended in recovery medium from Gibco (Cat. No 12648-010).

### **RECEIPT AND STORAGE**

Upon receipt confirm that adequate dry-ice is present, and the cells are frozen. Immediately transfer to -80°C storage. Cells should be stored at -80°C or at lower temperature and are stable as supplied until the expiry date shown. Cells should be diluted and plated immediately after thawing.

### **BACKGROUND**

Currently, the AAV product development and approval processes have multiple challenges that require the development of new tools (1). The *iLite*<sup>®</sup> AAV8 Packaging Assay Ready Cells together with the *iLite*<sup>®</sup> AAV Responsive Reporter Assay Ready Cells represent a functional, robust, and agile platform to facilitate the determination and titration of neutralizing antibodies against AAV vectors of serotype 8.

### **APPLICATION**

The *iLite*<sup>®</sup> AAV8 Packaging Assay Ready Cells together with the *iLite*<sup>®</sup> AAV Responsive Reporter Assay Ready Cells can be used as a transduction inhibition (TI) assay for the determination and titration of neutralizing antibodies against AAV vectors of serotype 8.

Application Notes for the following assays are available:

- Determination of NABs against AAV in human serum using *iLite*<sup>®</sup> AAV Packaging and *iLite*<sup>®</sup> AAV Responsive Reporter Assay Ready Cells (LABEL-DOC-00580)
- Titration of NABs against AAV in human serum using *iLite*<sup>®</sup> AAV Packaging and *iLite*<sup>®</sup> AAV Responsive Reporter Assay Ready Cells (LABEL-DOC-00599)







### **RELATED PRODUCTS**

<b>REF</b>	<b>Product name</b>
BM6100	<i>iLite</i> <sup>®</sup> AAV Responsive Reporter Assay Ready Cells
BM6002	<i>iLite</i> <sup>®</sup> AAV2 Packaging Assay Ready Cells
BM6005	<i>iLite</i> <sup>®</sup> AAV5 Packaging Assay Ready Cells
BM6006	<i>iLite</i> <sup>®</sup> AAV6 Packaging Assay Ready Cells
BM6009	<i>iLite</i> <sup>®</sup> AAV9 Packaging Assay Ready Cells

## REFERENCES

1. Gupta V, Lourenço SP, Hidalgo IJ. Development of Gene Therapy Vectors: Remaining Challenges. J Pharm Sci. 2020

## SYMBOLS ON LABEL

	Lot number		Temperature limitation
	Catalogue number		Biological risk
	Use by		Manufacturer

## PRECAUTIONS

For research use only. This product is intended for professional laboratory research use only. The data and results originating from using the product should not be used either in diagnostic procedures or in human therapeutic applications.

*iLite*<sup>®</sup> AAV8 Packaging Assay Ready Cells are a cell line of human origin classified as a Class 2 Genetically Modified Microorganism. They should be handled in accordance with EU directive (2009/41/EC) and disposed of in a licensed contained-use facility in accordance with these regulations. When used in accordance with the manufacturer's product specification, the requirements of EC Directive 2009/41/EC on the contained-use of genetically modified microorganisms are deemed to have been met.

Residues of chemicals and preparations generally considered as biohazardous waste should be inactivated prior to disposal by autoclaving or using bleach. All such materials should be disposed of in accordance with established safety procedures.

## PROPRIETARY INFORMATION

In accepting delivery of *iLite*<sup>®</sup> Assay Ready Cells the recipient agrees not to sub-culture these cells, attempt to sub-culture them or to give them to a third party, and only to use them directly in assays. *iLite*<sup>®</sup> cell-based products are covered by patents which is the property of Svar Life Science AB and any attempt to reproduce the delivered *iLite*<sup>®</sup> Assay Ready Cells is an infringement of these patents.

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES

