

## HyClone™ Peak Expression Medium

## Instructions for Use

## **Product description**

HyClone<sup>TM</sup> peak expression medium (the product) is intended to be used for both growth and transfection of human embryonic kidney cells (HEK293) for adeno-associated virus (AAV) production.

Peak expression medium is an animal-derived component-free (ADCF), hydrolysate-free, and serum-free cell culture medium. This product contains a stable form of glutamine and poloxamer 188, and supports for high-yield virus production and increased cell culture efficiency.

This product is available as ready-to-use liquid medium, or as powder medium.

## Safety

For use and handling of the product, refer to the Safety Data Sheet on *cytiva.com*.

## **Storage**

Liquid and powder medium should be stored at  $2^{\circ}$ C to  $8^{\circ}$ C in a dry environment, protected from light. In addition, powder medium should be stored in a tightly sealed container to protect from moisture.

#### **Shelf life**

Liquid medium is stable up to 12 months and powder medium is stable up to 24 months from the date of manufacture, when stored at 2°C to 8°C and protected from light.

Note: Refer to the product label for the expiry date.

## **Disposal**

Dispose contents and container in accordance with all local and national regulations.

### **Specifications and components**

The specifications of this product are as follows:

- ADCF formulation
- Hydrolysate-free
- Serum-free

The components for the liquid and powder medium are listed below.

Components	Liquid medium	Powder medium
Alanyl-glutamine (4 mM)	+	+
Poloxamer 188 (0.08%)	+	+
Sodium bicarbonate	+	-
Phenol red	-	-

# Required materials for hydration of powder medium

The following equipment and materials, not provided with the product, are required to hydrate the powder medium.

- Mixing vessel
- Stirrer

**Note:** A magnetic stirrer can be used for small-scale hydration up to 5 L. An overhead or bottom-mounted impeller is recommended for larger volumes.

- Calibrated pH meter
- Calibrated osmometer
- Cell culture-grade water
- Sodium bicarbonate
- 0.2 µm filter
- Container for hydrated medium

## **Hydration of powder medium**

Follow the steps below to hydrate the powder medium.

Ste	n	Acti	ion
<b>-</b>	~	700	

- Fill a clean mixing vessel with 90% of the final volume with cell culture-grade water at ambient temperature (20°C to 25°C). For example, start with 900 mL of water for a preparation of 1 L medium. Start stirring.
- Slowly add 25.17 g/L powder to the vessel to avoid clumping. Mix for 20 minutes until dissolved.
- 3 Slowly add 2.50 g/L sodium bicarbonate. Mix for 10 minutes until all components are completely dissolved.

Result.

The medium should be a pale yellow and particulate-free solution. Prior to pH adjustment, the medium can be slightly hazy.

cytiva.com 29717925 AB

#### Step Action

4 Adjust the pH to 7.2 (acceptable range pH 7.0 to 7.4) by dropwise addition of 5 N NaOH or HCI.



#### **NOTICE**

Use caution when adjusting pH. Over adjusting can cause the osmolality to be out of specification.

- 5 Adjust the medium to the final volume with cell culture-grade water at ambient temperature (20°C to 25°C).
- 6 Measure and record the final pH and osmolality.

Final values must be:

- pH 7.0 to 7.4
- Osmolality 300 to 340 mOsm/kg
- 7 Sterile filter the medium immediately after hydration using a 0.2 µm filter.
- 8 Store the hydrated medium in a container, protected from light at 2°C to 8°C until use.

#### General culture recommendations

For optimal culture performance, cultures should be incubated at 37°C in a 5% CO<sub>2</sub> environment.

#### **Cell adaptation**

Adaptation of HEK293 cells from another medium to this product can be achieved by two methods:

- Direct adaptation
- Gradual adaptation

Follow the recommendations below for direct or gradual adaptation of HEK293 cells.

### **Direct adaptation**

- Seed suspension HEK293 cells from the current medium to this product at a cell density of at least 3×10<sup>5</sup> viable cells/mL. Increasing the seeding density 2 to 3 times (i.e., 6×10<sup>5</sup> viable cells/mL or 9×10<sup>5</sup> viable cells/mL) can be beneficial.
- Passage cells every 3 to 4 days.
- Passage cells 3 to 4 times in this product to make sure that the adaptation is complete. The cells should have at least 90% viability after 3 to 4 passages.
- If the cell viability does not reach 90% after 3 to 4 passages, consider switching to a gradual adaptation approach.
- If the cell density is low at scheduled passage intervals, consider waiting another 24 to 36 hours to passage.
- Make sure that cells are completely adapted to this product before performing screening and/or benchmark studies.

#### **Gradual adaptation**

 Mix the current medium and this product at a stepwise ratio of 90:10, 75:25, 50:50, 25:75, and 0:100.

**Tip:** Customized combinations can also be beneficial.

- Seed suspension HEK293 cells from the current medium to this product at a cell density of at least 3×10<sup>5</sup> viable cells/mL.
  Increasing the seeding density 2 to 3 times can be beneficial at later ratios (50:50, 25:75, etc.).
- Make sure that cells are healthy and reach normal densities prior to moving to the next step of adaptation.
- Cells can be spun down to remove spent media prior to passage into fresh medium (optional).

**Note:** Conditioned media carried over during passaging can be beneficial for certain cell lines.

- Passage cells 3 to 4 times in this product to make sure that the adaptation is complete. The cells should have at least 90% viability after 3 to 4 passages.
- Make sure that cells are completely adapted to this product before performing screening and/or benchmark studies.

#### Cryopreservation

Adapted cells can be cryopreserved in this product with 10% Dimethyl Sulfoxide (DMSO). The recommended cell density for freezing is a minimum of  $1.0 \times 10^7$  viable cells/mL.

### **Quality control testing**

Quality control test specifications are listed in the table below.

Parameter	Specification	
Appearance	Powder medium: off-white powder	
	Liquid medium: pale yellow solution	
pH (at 20°C to 25°C)	7.0 to 7.4 (with NaHCO <sub>3</sub> )	
Osmolality	$300  \mathrm{to}  340  \mathrm{mOsm/kg}  \mathrm{H}_2\mathrm{O}$	
Sterility	No growth (bacteria or fungi)	
Endotoxin	< 1.0 EU/mL	
HPLC amino acid profile	Complies (80% to 120%) recovery	

Refer to Certificate of analysis for actual results.

## **Ordering information**

Product	Pack size	Product code
HyClone peak expression powder medium	5 L HDPE bottle	SH31192.01
	10 L HDPE bottle	SH31192.02
	50 L HDPE bottle	SH31192.03
	100 L poly bag/pail	SH31192.04
	500 L poly bag/pail	SH31192.05
	1000 L poly bag/drum	SH31192.06
HyClone peak expression liquid medium	1 L bottle	SH31193.02
	1 L bag	SH31193.08
	5 L bag	SH31193.09
	100 L bag	SH31193.10
	200 L bag	SH31193.11
	10 L bag	SH31193.12
	20 L bag	SH31193.13
	50 L bag	SH31193.14

#### **Technical support**

Contact your local Cytiva representative to learn more about the services we offer. To find a certificate or SDS for a specific product, visit cytiva.com/certificates.





## $\textbf{Give feedback} \, \textbf{on this document} \,$

Visit cytiva.com/techdocfeedback or scan the QR code.



## cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

 $Hy Clone is a trademark of Global Life Sciences Solutions \, USA \, LLC \, or \, an \, affiliate \, doing \, business \, as \, Cytiva.$ 

Any other third-party trademarks are the property of their respective owners.

© 2022–2023 Cytiva

For local office contact information, visit cytiva.com/contact 29717925 AB V:5 01/2023

