

HyClone™ glycan supplement

Product Information

Product description

HyClone™ glycan supplement is a chemically defined and animal-derived component-free (ADCF) cell culture supplement. It is designed to modulate galactosylation profiles of therapeutic proteins during fed-batch culture of Chinese hamster ovary (CHO) cell lines.

The product is supplied as a ready-to-use liquid supplement and is available in pack sizes suitable for both small-scale and large-scale bioprocessing applications.

For optimal performance, the product is recommended for use in combination with the HyClone cell culture platform, including the following:

- HyClone ActiPro™ medium
- HyClone Cell Boost™ 7a supplement
- HyClone Cell Boost 7b supplement

The product is hydrolysate-free and serum-free and is designed to support efficient cell culture performance and protein production.

Compatibility

The product has been validated for use in CHO cell lines. Performance and compatibility in other cell lines have not been evaluated.

Safety

Use aseptic techniques when handling the product. For the details regarding use and handling of the product, refer to the *Safety Data Sheet* (SDS) on [cytiva.com](https://www.cytiva.com).

Storage

Store the product at 2°C to 8°C in a tightly sealed container.

Shelf life

The product is stable up to 6 months from the date of manufacture, when stored under the conditions specified above.

Note: For the expiry date, refer to the product label.

Disposal

Dispose of the product container and its contents in accordance with all local and national regulations.

General cell culture recommendations

For optimal cell culture performance, incubate CHO cells under the conditions recommended for the specific cell line. For additional guidance on CHO cell culture, refer to the data files for HyClone ActiPro medium, HyClone Cell Boost 7a supplement, and HyClone Cell Boost 7b supplement, available on [cytiva.com](https://www.cytiva.com).

Supplementation recommendations

Start supplementation on day 3 at 0.5% (v/v). Continue supplementing the culture at the same concentration every other day according to the schedule in the table below. The schedule serves as a baseline and is adjustable to meet specific process requirements and glycosylation targets.

Days	Supplementation
3, 5, 7, 9, and 11	0.5% (v/v)

Harvest recommendation

It is recommended to harvest the cells before the cell culture drops below 70% viability.

Quality control testing

Quality control test specifications are listed in the table below.

Parameter	Specification
Appearance	Clear, colorless solution
Osmolality	285–325 mOsm/kg H ₂ O (1:4 dilution)
Sterility	No growth (bacteria or fungi)
Endotoxin	< 1.0 EU/mL

For actual results, refer to the *Certificate of Analysis*, available on [cytiva.com/certificates](https://www.cytiva.com/certificates).

Ordering information

Product	Pack size	Product code
HyClone glycan supplement	100 mL bottle*	SH31210.01
	1000 mL bottle	SH31210.02

* Bottles are made of polyethylene terephthalate (PET).

Related product	Pack size	Product code
HyClone ActiPro medium, liquid	500 mL bottle	SH31039.01
	1000 mL bottle	SH31039.02
	1 L bottle	SH31039.03
HyClone ActiPro medium, dry powder	5 L bottle	SH31037.01
	10 L bottle	SH31037.02
	25 L bottle	SH31037.05
HyClone Cell Boost 7a, liquid	500 mL bottle	SH31119.01
	1000 mL bottle	SH31119.02
	1 L bag	SH31119.03
	5 L bag	SH31119.04
	10 L bag	SH31119.05
	20 L bag	SH31119.06
	50 L bag	SH31119.07
	100 L bag	SH31119.08
	200 L bag	SH31119.09
HyCloneCell Boost 7a, dry powder	1 L bottle	SH31026.07
	5 L bottle	SH31026.01
	10 L bottle	SH31026.02
	25 L bottle	SH31026.03

Related product	Pack size	Product code
HyClone Cell Boost 7b, liquid	100 mL bottle	SH31120.01
	500 mL bottle	SH31120.02
	1000 mL bottle	SH31120.03
	1 L bag	SH31120.04
	5 L bag	SH31120.05
	10 L bag	SH31120.06
	20 L bag	SH31120.07
	50 L bag	SH31120.08
	100 L bag	SH31120.09
	200 L bag	SH31120.10
HyClone Cell Boost 7b, dry powder	0.5 L bottle	SH31027.01
	1 L bottle	SH31027.07
	5 L bottle	SH31027.02
	10 L bottle	SH31027.04

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](https://www.cytiva.com/certificates).



Give feedback on this document

Visit [cytiva.com/techdocfeedback](https://www.cytiva.com/techdocfeedback) or scan the QR code.



cytiva.com

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

ActiPro, Cell Boost, HyClone, and HyClone Cellboost are trademarks 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.

© 2026 Cytiva

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

30029986 AA V:2 04/2026

