

Certificate of Analysis

CloneMedia CHO Growth A, without L-Gln K8840

Batch #	99901150703
----------------	--------------------

Full Description	Semi-solid media for the growth of suspension-adapted CHO, DG44 and DUXB11 cells. Contains methylcellulose, recombinant proteins and other supplements
Appearance	Viscous liquid frozen for storage and transport
Components	No material of animal origin is used in the manufacture of this product
Volume	90ml for 100ml of media
Sterility	Pass. Note: USP <71> Direct Inoculation method used.
Preservatives/Stabilisers	None
Endotoxin Levels	<0.05 EU/ml endotoxin unit (EU) per ml as determined by LAL method
Functional Growth	Supports growth of in house cell line
Cell Line Used	CHO-S Cell Line (Life Technologies, Cat # 11619-012)
Expiry Date	Up to 6 months from receipt when stored at -20°C Once defrosted store at 2°C - 8°C and use within 1 week

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

For research use only. Store at -20°C

Signed: *Kieu-Khiem Nguyen*

Date: 04 SEP 2015

Production Biologist