

# Certificate of Analysis

## CloneMedia CHO Growth A, without L-Gln K8840

<b>Batch #</b>	<b>99901171204</b>
----------------	--------------------

<b>Full Description</b>	Semi-solid media for the growth of suspension-adapted CHO, DG44 and DUXB11 cells. Contains methylcellulose, recombinant proteins and other supplements
<b>Appearance</b>	Viscous liquid frozen for storage and transport
<b>Components</b>	No material of animal origin is used in the manufacture of this product
<b>Volume</b>	90ml for 100ml of media
<b>Sterility</b>	Pass. Note: USP <71> Direct Inoculation method used.
<b>Preservatives/Stabilisers</b>	None
<b>Endotoxin Levels</b>	<.05 EU/ml endotoxin unit (EU) per ml as determined by LAL method
<b>Functional Growth</b>	Supports growth of in house cell line
<b>Cell Line Used</b>	CHO-S Cell Line (Life Technologies, Cat # 11619-012)
<b>Expiry Date</b>	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 ISO9001:2015 Quality Assurance guidelines

**For research use only. Store at -20°C**

Signed: *Kieu-Khiem Nguyen*

Date: 21-FEB-2018

**Reagent Biologist**