

Certificate of Analysis

CloneMedia CHO Growth A, without L-Gln K8840

| | |
|----------------|--------------------|
| Batch # | 99901190801 |
|----------------|--------------------|

| | |
|----------------------------------|--|
| 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 | <.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 ISO9001:2015 Quality Assurance guidelines

For research use only. Store at -20°C

Signed: *Kieu-Khiem Nguyen*

Date: 16-OCT-2019

Reagent Biologist