

Certificate of Analysis

CloneXL

K8521

| | |
|----------------|-----------------|
| Batch # | CXL0715A |
|----------------|-----------------|

| | |
|----------------------------------|---|
| Full Description | Animal-free, chemically-defined supplement for CHO-S |
| Appearance | Clear liquid, frozen for storage and transport |
| Components | No material of animal origin is used in the manufacture of this product |
| Volume | 2ml for 100ml of media |
| Sterility | No bacterial or fungal growth after 72 hours incubation |
| Preservatives/Stabilisers | None |
| Endotoxin Levels | <0.25 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 Invitrogen Cat # 11619-012 |
| Expiry Date | 12 months from receipt if stored at -20°C Once defrosted use immediately |

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

For research use only Store at -20°C

Signed: *Kieu-Khiem Nguyen*

Date: 09 SEP 2015

Production Biologist