

# Certificate of Analysis

## CloneXL

## K8521

|                |                 |
|----------------|-----------------|
| <b>Batch #</b> | <b>CXL1014A</b> |
|----------------|-----------------|

|                                  |   |
|----------------------------------|---|
| <b>Full Description</b>          | Animal-free, chemically-defined supplement for CHO-S                        |
| <b>Appearance</b>                | Clear 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>                    | 2ml for 100ml of media  |
| <b>Sterility</b>                 | No bacterial or fungal growth after 72 hours incubation                     |
| <b>Preservatives/Stabilisers</b> | None  |
| <b>Endotoxin Levels</b>          | <0.100 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<br>Invitrogen Cat # 11619-012                               |
| <b>Date of manufacture</b>       | 03-SEP-2014   |
| <b>Expiry Date</b>               | 12 months from receipt if stored at -20°C<br>Once defrosted use immediately |

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

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

Signed: *Christopher Lee*

Date: 22-OCT-2014

**Production Biologist**