

# Certificate of Analysis

## CloneXL K8521

<b>Batch #</b>	<b>CXL1020A</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>	<0.5 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 Invitrogen Cat # 11619-012
<b>Expiry Date</b>	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: 25-MAR-2021

**Production Biologist**