

Certificate of Analysis

CloneXL

K8521

Batch #	CXL0113A
----------------	----------

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	<0.100 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
Date of manufacture	04 January 2013
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:

Date: 7 February 2013

Production Supervisor