

# Certificate of Analysis

## CloneMedia-CHO

### K8710

<b>Batch #</b>	CMC0612A
----------------	----------

<b>Full Description</b>	Semi-solid media for the growth of suspension-adapted CHO cells, contains methylcellulose, stable glutamine (8mM) and supplements
<b>Appearance</b>	Viscous 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>	90ml 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 Invitrogen Cat # 11619-012
<b>Date of manufacture</b>	10 May 2012
<b>Expiry Date</b>	Up to 12 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 ISO9001:2008 Quality Assurance guidelines

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



Signed:

Date: 26 July 2012

**Production Supervisor**