

Certificate of Analysis

CloneMedia-CHO-G DHFR K8742

Batch #	CMD0914A
----------------	-----------------

Full Description	Semi-solid media for the growth of suspension-adapted CHO, DG44 and DUXB11 cells under GS selection, contains methylcellulose, recombinant proteins and supplements
Appearance	Viscous liquid frozen for storage and transport
Components	No material of animal origin is used in the manufacture of this product
Volume	90ml for 100ml of media
Sterility	No bacterial or fungal growth after 72 hours incubation
Preservatives/Stabilisers	None
Endotoxin Levels	< 0.3 EU/ml endotoxin unit (EU) per ml as determined by LAL method
Functional Growth	Supports growth of in house cell line (L-Glutamine supplanted)
Cell Line Used	CHO-S Cell Line Invitrogen Cat # 11619-012
Date of manufacture	17-Sep-2014
Expiry Date	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: 06-Oct-2014



Production Biologist