

Certificate of Analysis

CloneMedia-CHOK1SV

K8725

Batch #	CMV0114A
----------------	-----------------

Full Description	Glutamine free semi-solid media for the growth of suspension-adapted CHOK1SV cells. Contains methylcellulose, recombinant proteins and other 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
Endotoxin Levels	0.441 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	Sep 20 th , 2013
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: 11-Mar-2014

Christopher Lee

Production Biologist