

Certificate of Analysis K8510 CloneMatrix

Batch #	MCW0611A
Full Description	Methylcellulose in DI Water
Appearance	Viscous liquid frozen for storage and transport
Components	No material of animal origin is used in the manufacture of this product
Volume	40ml for 100ml of media
Sterility	No bacterial or fungal growth after 72 hours incubation
Preservatives/Stabilizers	None
Endotoxin levels	0.1 EU/ml endotoxin unit (EU) per ml as determined by LAL method
Functional Growth	Supports growth of in house cell line
Cell line used	Hybridoma Cell Line ECACC Cat. # 90112228
Date of manufacture	June 2011
Expiry Date	Up to three years from receipt if stored at -20°C Once defrosted store at 2-8°C for up to 6 months.

This product is manufactured to ISO13485:2003/CMDCAS certification.
Manufactured in an FDA registered facility compliant with QSR
For Research use only. Store at -20°C

Signed: *R Mitchell*

Date: 28/06/11

Production Biologist

