

Certificate of Analysis

CloneMatrix

K8510

Batch #	MCW0417A
----------------	-----------------

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	0.2µm sterile filtered
Preservatives/Stabilisers	None
Endotoxin Levels	<0.50 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
Expiry Date	Up to three years from receipt if stored at -20°C Once defrosted store at 2°C – 8°C for up to 6 months

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

For research use only Store at -20°C

Signed: Kieu Khiem Nguyen

Date: 23-JUN-2017

Production Biologist