

# Certificate of Analysis

## CloneMatrix

### K8510

<b>Batch #</b>	<b>MCW0619A</b>
----------------	-----------------

<b>Full Description</b>	Methylcellulose in DI Water
<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>	40ml for 100ml of media
<b>Sterility</b>	0.2µm sterile filtered
<b>Preservatives/Stabilisers</b>	None
<b>Endotoxin Levels</b>	<0.5 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>	Hybridoma Cell Line ECACC Cat # 90112228
<b>Expiry Date</b>	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: 01-JUL-2019*

**Production Biologist**