

Certificate of Analysis

CloneMatrix

K8510

| | |
|----------------|-----------------|
| Batch # | MCW1117A |
|----------------|-----------------|

| | |
|----------------------------------|---|
| 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 | <0.5 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: 22-NOV-2017

Production Biologist