

# Certificate of Analysis

## CloneMatrix K8510MD

|                |                |
|----------------|----------------|
| <b>Batch #</b> | <b>1495862</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<br>ECACC Cat # 90112228   |
| <b>Expiry Date</b>               | Up to three years from receipt if stored at -20°C<br>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: *Erwin Ramirez*

Date: 31-May-2018

**Production Biologist**