

Certificate of Analysis

CloneMedia

K8610

Batch #	CMA1215A
----------------	-----------------

Full Description	Semi-solid media for the growth of hybridoma/myeloma
Appearance	Viscous liquid frozen for storage and transport
Components	Contains methylcellulose, FBS, stable L-Glutamine and additional supplements
Volume	90ml for 100ml of media
Sterility	No bacterial or fungal growth after 72 hours incubation
Functional Growth	Supports growth of in house cell line
Cell Line Used	Primary mouse spleen cells fused with P3x63Ag8.653 myeloma cells (ATCC Cat # CRL1580)
Endotoxin Levels	< 2.5 EU/ml endotoxin unit (EU) per ml
Expiry Date	Up to 12 months from receipt when stored at -20°C Once defrosted store at 2°C – 8°C and use within 1 week

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

For research use only Store at -20°C

Signed: *Kieu-Khiem Nguyen*

Date: 26-JAN-2016

Production Biologist