

Certificate of Analysis

CloneMedia Hybridoma Semi-Solid Selection and Cloning Medium (with HAT)

K8865

Lot #	17H83484
--------------	-----------------

Full Description	Semi-solid methylcellulose-based medium for selecting and cloning of hybridoma clones.
Appearance	Viscous, red liquid frozen for storage and transport
Components	Contains methylcellulose, Dulbecco's Modified Eagle's Medium, FBS, Hypoxanthine, aminopterin, thymidine (HAT) and additional supplements
Volume	90 mL for 100 mL of medium
Sterility	Pass (No growth after 14 days)
Preservatives/Stabilisers	None
Functional Growth	Pass (Cell culture performance >80% compared to standard)
Cell Line Used	Hybridoma cell line
Expiry Date	6 months from date of receipt

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

For research use only. Store at -20°C

Signed:

Production



Date:

9/29/2017