

# Certificate of Analysis

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

### K8865

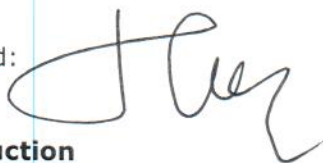
<b>Lot #</b>	<b>1000014060</b>
--------------	-------------------

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

1/02/2020