

Certificate of Analysis

XP Media Hybridoma Fusion Recovery Medium K8864

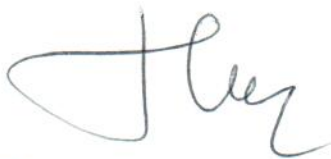
Lot #	17M86921
--------------	-----------------

Full Description	Medium formulated to promote hybridoma viability following the fusion process, but prior to clone selection
Appearance	Clear, red liquid frozen for storage and transport
Components	Contains Dulbecco's Modified Eagle's Medium, FBS and additional supplements
Volume	100 mL
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:



Date:

1/22/2018

Production