

Certificate of Analysis

XP Media Hybridoma Fusion Medium

K8863

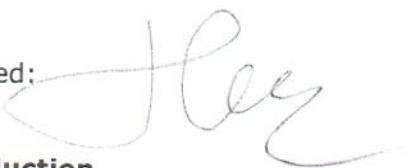
Lot #	17D78702
--------------	-----------------

Full Description	Medium for washing lymphocytes and parental myeloma cells to remove serum prior to fusion and for use during the fusion process
Appearance	Clear, red liquid frozen for storage and transport
Components	Contains Dulbecco's Modified Eagle's Medium and additional supplements
Volume	500 mL
Sterility	Pass (No growth after 14 days)
Preservatives/Stabilisers	None
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:

11/27/2017

Production