

Certificate of Analysis

XP Media Pre-Fusion Myeloma Growth Medium and Hybridoma Expansion Medium (without HAT)

K8862

Lot #	17B78541
--------------	-----------------

Full Description	Complete medium for the growth of myeloma cells prior to fusion and to support the expansion of hybridoma clones
Appearance	Clear, red liquid frozen for storage and transport
Components	Contains Dulbecco's Modified Eagle's Medium, FBS 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: 3/27/2017

Production