

Certificate of Analysis

XP Media Hybridoma Growth Medium (with HT)

K8866

Lot #	16E71279
-------	----------

Full Description	Rich, complete medium optimized for hybridoma expansion following clone selection and colony picking
Appearance	Clear, red liquid frozen for storage and transport
Components	Contains Dulbecco's Modified Eagle's Medium, FBS, hypoxanthine, thymidine (HT) and additional supplements
Volume	500 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:

6/13/16

Production

