

Certificate of Analysis

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

K8862

| | |
|--------------|-----------------|
| Lot # | 18E91191 |
|--------------|-----------------|

| | |
|----------------------------------|--|
| 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:

9/24/2019

Production