

# Certificate of Analysis

## XP Media Hybridoma Fusion Medium

### K8863

|       |          |
|-------|----------|
| Lot # | 16E71272 |
|-------|----------|

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



**Production**

Julie Nguyen

Date:

6/13/16