

# Certificate of Analysis

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


### K8862

<b>Lot #</b>	<b>1000003101</b>
--------------	-------------------

<b>Full Description</b>	Complete medium for the growth of myeloma cells prior to fusion and to support the expansion of hybridoma clones
<b>Appearance</b>	Clear, red liquid frozen for storage and transport
<b>Components</b>	Contains Dulbecco's Modified Eagle's Medium, FBS and additional supplements
<b>Volume</b>	500 mL
<b>Sterility</b>	Pass (No growth after 14 days)
<b>Preservatives/Stabilisers</b>	None
<b>Expiry Date</b>	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: 10/21/2019

**Production**