

# Certificate of Analysis

## XP Media Hybridoma Fusion Recovery Medium K8864

<b>Lot #</b>	<b>18L96979</b>
--------------	-----------------

<b>Full Description</b>	Medium formulated to promote hybridoma viability following the fusion process, but prior to clone selection
<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>	100 mL
<b>Sterility</b>	Pass (No growth after 14 days)
<b>Preservatives/Stabilisers</b>	None
<b>Functional Growth</b>	Pass (Cell culture performance >80% compared to standard)
<b>Cell Line Used</b>	Hybridoma cell line
<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:

12/20/2018

**Production**