

# Certificate of Analysis

## XP Media CHO Growth A without L-Gln K8860

<b>Batch #</b>	<b>99914210101</b>
----------------	--------------------

<b>Full Description</b>	XP Media CHO Growth A without L-glutamine
<b>Appearance</b>	Light pink solution
<b>Components</b>	No material of animal origin is used in the manufacture of this product
<b>Volume</b>	1000 ml of media
<b>Sterility</b>	Pass. Note: USP <71> Direct Inoculation method used.
<b>Preservatives/Stabilisers</b>	None
<b>Endotoxin Levels</b>	<.05 EU/ml endotoxin unit (EU) per ml as determined by LAL method
<b>Functional Growth</b>	Supports growth of in house cell line

This product has been manufactured to ISO9001:2015 Quality Assurance guidelines

**For research use only. Store at 2-8°C**

Signed: *Kieu Khiem Nguyen*

Date: 22-FEB- 2021

**Reagent Biologist**