

# Certificate of Analysis

## XP Media-CHO-G

### K8760

<b>Batch #</b>	<b>CMF1212A</b>
----------------	-----------------

<b>Full Description</b>	Liquid media for the growth of suspension-adapted CHO cells under GS selection, animal-free, chemically-defined
<b>Appearance</b>	Clear yellow solution frozen for storage and transport
<b>Components</b>	No material of animal origin is used in the manufacture of this product
<b>Volume</b>	500ml media
<b>Sterility</b>	No growth of aerobic or anaerobic bacteria or fungi detected
<b>Preservatives/Stabilisers</b>	None
<b>pH</b>	7.3
<b>Osmolarity</b>	312 mOsmol/kg
<b>Endotoxin Levels</b>	0.16 endotoxin unit (EU) per ml
<b>Functional Growth</b>	Supports growth of in house cell line (L-Glutamine supplemented)
<b>Cell Line Used</b>	CHO-S Cell Line Invitrogen Cat # 11619-012
<b>Date of manufacture</b>	17-Jul-2012
<b>Expiry Date</b>	Up to 24 months from receipt when stored at -20°C Once defrosted store at 2°C – 8°C and use within 6 months

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

**For research use only    Store at -20°C**

Signed:

Date: 20-Dec-2012

*Christopher Lee*

**Production Chemist**