

Certificate of Analysis

XP Media-CHO

K8750

Batch #	CME1212A
----------------	-----------------

Full Description	Liquid media for the growth of suspension-adapted CHO cells, animal-free, chemically-defined, contains stable glutamine (8mM)
Appearance	Clear yellow solution frozen for storage and transport
Components	No material of animal origin is used in the manufacture of this product
Volume	500ml media
Sterility	No growth of aerobic or anaerobic bacteria or fungi detected
Preservatives/Stabilisers	None
pH	7.2
Osmolarity	322 mOsmol/kg
Endotoxin Levels	0.12 endotoxin unit (EU) per ml
Functional Growth	Supports growth of in house cell line
Cell Line Used	CHO-S Cell Line Invitrogen Cat # 11619-012
Date of manufacture	20-Jul-2012
Expiry Date	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 Chemsit