

Certificate of Analysis

XP Media CHO Growth A without L-Gln K8860

Batch #	99914151001
----------------	--------------------

Full Description	XP Media CHO Growth A without L-glutamine
Appearance	Light pink solution
Components	No material of animal origin is used in the manufacture of this product
Volume	1000 ml of media
Sterility	Pass. Note: USP <71> Direct Inoculation method used.
Preservatives/Stabilisers	None
Endotoxin Levels	<.03 EU/ml endotoxin unit (EU) per ml as determined by LAL method
Functional Growth	Supports growth of in house cell line
Expiry Date	Stated on the bottle when stored at 2-8°C

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

For research use only. Store at 2-8°C

Signed: *Kieu-Khiem Nguyen*

Date: 12 MAY 2016

Production Biologist