

Certificate of Analysis

XP Media CHO Growth A without L-Gln K8860

Batch #	99914181201
----------------	--------------------

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	<.05 EU/ml endotoxin unit (EU) per ml as determined by LAL method
Functional Growth	Supports growth of in house cell line
Expiry Date	Up to 6 months from receipt when stored at 2-8°C

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: 29 JAN 2019

Production Biologist