

Certificate of Analysis

XP Media

K8650

Batch #	CMJ0812A
----------------	----------

Full Description	Liquid media for hybridoma and myeloma cells
Appearance	Clear red solution frozen for storage and transport
Components	FBS (PAA Product: A11-204 Lot: A20410-2463 L-Glutamine (2mM) and supplements
Volume	500ml media
Sterility	No growth of aerobic or anaerobic bacteria or fungi detected
pH	7.1
Osmolarity	365 mOsmol/kg
Endotoxin Levels	0.83 endotoxin unit (EU) per ml
Functional Growth	Supports growth of in house cell line
Cell Line Used	Hybridoma Cell Line ECACC Cat # 90112228
Date of manufacture	30 April 2012
Expiry Date	Up to 12 months from receipt when stored at -20°C Once defrosted store at 2°C – 8°C and use within 3 months

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

For research use only Store at -20°C



Signed:

Date: 21 August 2012

Production Supervisor