
Certificate of AnalysisK8690
XP Media-HEK 293**Genetix****Batch #** CME0108/71260

Full Description	Liquid media for the growth of suspension-adapted HEK cells.
Appearance	Clear solution frozen for storage and transport
Components	Media may contain discrete proteins or bulk protein fractions.
Volume	500ml media
Sterility	No growth of aerobic bacteria, anaerobic bacteria or fungi detected
Preservatives/Stabilizers	BSA (PAA K41-001, Lot # K00107-0714)
pH	7.1 – 7.5
Osmolality	267 Omol/kg
Endotoxin levels	<0.05 endotoxin unit (EU) per ml
Functional Growth	Supports growth of in house cell line
Cell line used	HEK 293 - F cell Line (Invitrogen # R790-07)
Date of manufacture	28 th November 2007

This product is manufactured in an FDA registered, cGMP compliant facility

For Research use only. Store at -20°C.

Signed _____ Date _____

Reagents Laboratory Manager