

Certificate of Analysis

CloneMedia-HEK

K8685

| | |
|----------------|-----------------|
| Batch # | CMI0313A |
|----------------|-----------------|

| | |
|----------------------------------|--------------------------------------------------------------------------------------------------------------|
| Full Description | Semi-solid media for the growth of suspension-adapted HEK cells |
| Appearance | Viscous liquid frozen for storage and transport |
| Components | No material of animal origin is used in the manufacture of this product |
| Volume | 90ml for 100ml of media |
| Sterility | No bacterial or fungal growth after 72 hours incubation |
| Preservatives/Stabilisers | None |
| Endotoxin Levels | <0.3 EU/ml endotoxin unit (EU) per ml as determined by LAL method |
| Functional Growth | Supports growth of in house cell line |
| Cell Line Used | Freestyle HEK 293 Cell Line Freestyle Cat # R790-07 |
| Date of manufacture | 22 March 2013 |
| Expiry Date | Up to 12 months from receipt when stored at -20°C Once defrosted store at 2°C - 8°C and use within 1 week |

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

For research use only Store at -20°C



Signed:

Date: 11 April 2013

Production Supervisor