

# Certificate of Analysis

## CloneMedia-HEK

### K8685

<b>Batch #</b>	<b>CM10620A</b>
----------------	-----------------

<b>Full Description</b>	Semi-solid media for the growth of suspension-adapted HEK cells
<b>Appearance</b>	Viscous liquid frozen for storage and transport
<b>Components</b>	No material of animal origin is used in the manufacture of this product
<b>Volume</b>	90ml for 100ml of media
<b>Sterility</b>	No bacterial or fungal growth after 72 hours incubation
<b>Preservatives/Stabilisers</b>	None
<b>Endotoxin Levels</b>	0.1772 EU/ml endotoxin unit (EU) per ml as determined by LAL method
<b>Functional Growth</b>	Supports growth of inhouse cell line
<b>Cell Line Used</b>	Freestyle HEK 293 Cell Line Freestyle Cat # R790-07
<b>Expiry Date</b>	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: *Kieu-Khiem Nguyen* Date: 25-JUN-2020

**Production Biologist**