

# Certificate of Analysis

## CloneMedia-PER.C6

### K8775

<b>Batch #</b>	<b>CMP0514A</b>
----------------	-----------------

<b>Full Description</b>	Semi-solid media for the growth of PER.C6® cells. Includes stable glutamine (4mM), recombinant proteins and supplements
<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>	< 0.3 EU/ml endotoxin unit (EU) per ml as determined by LAL method
<b>Functional Growth</b>	Supports growth of in house cell line
<b>Cell Line Used</b>	PER.C6 Cell Line (PER.C6® is a registered trademark of Crucell N.V)
<b>Date of manufacture</b>	14-May-2014
<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:

Date: 10-Jun-14

*Christopher Lee*

**Production Biologist**