

Certificate of Analysis

Hybridoma Polyethylene Glycol for Cell Fusion

K8868

Lot #	19E102728
--------------	------------------

Full Description	Solution used for the fusion of mouse splenocytes and parental myeloma cells to generate hybridomas
Appearance	Clear, pink liquid
Components	Contains Dulbecco's Modified Eagle's Medium and polyethylene glycol
Volume	1.5 mL
Sterility	Pass (No growth after 14 days)
Preservatives/Stabilisers	None
Expiry Date	6 months from date of receipt

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

For research use only. Store at 2-8°C

Signed:



Date:

9/30/19

Production