

Certificate of Analysis

Adherent Cell Picking Reagents Test Kit

Types A - D

K8005

Batch #	MD168845
----------------	-----------------

Full Description	Non-enzymatic buffered solution for picking of adherent mammalian cells
Appearance	Clear liquid
Components	No material of animal origin is used in the manufacture of this product
Volume	4 x 250ml
Buffer	PBS
Preservatives/Stabilisers	None
Sterility	Autoclaved at 121°C for 15 minutes
Date of manufacture	20-Mar-2013

This product has been manufactured to ISO9001:2008 certification.

For research use only Store at 2°C - 8°C

Christopher Lee

Signed:

Date: 20-Mar-2013

Production Biologist