

Certificate of Analysis

Adherent Cell Picking Reagent

Type B

K8020

Batch #	0813A
----------------	-------

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	1 x 500ml
Buffer	PBS
Preservatives/Stabilisers	None
Sterility	Autoclaved at 121°C for 15 minutes
Date of manufacture	29 August 2013

This product has been manufactured to ISO9001:2008 certification.

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



Signed:

Production Supervisor

Date: 29 August 2013