

Certificate of Analysis

Adherent Cell Picking Reagent

Type A

K8010

Batch #	MD158862
----------------	----------

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	09-APR-2012

This product has been manufactured to ISO9001:2008 certification.

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

Signed: 

Date: 09-APR-2012

Production Chemist