

Certificate of Analysis

CloneDetect

Human IgG (H+L) Specific, 550 Labelled K8210

Batch #	CDC1212A
----------------	-----------------

Full Description	Sheep anti-human IgG, polyclonal, F(ab)2, Dylight™ 550 conjugated. Sterile, azide free
Origin of Host Species	Sheep Serum
Immunogen	Human IgG
Format	F(ab)2 Conjugated to Dylight™ 550 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
Specificity	Human IgG
Method of Purification	Antigen affinity chromatography followed by pepsin digestion to obtain F(ab)2 fragments
Appearance	Clear pink solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	3.8
Buffer	Phosphate Buffered Saline pH 7.2
Preservatives/Stabilisers	None
Sterility	0.2µm sterile filtered

Date of manufacture	12-Decmeber-2012
Expiry Date	12 months from receipt of product

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

For research use only Store at 2 - 8°C

Signed:



Date: 26 December 2012

Production Supervisor