

# Certificate of Analysis

## CloneDetect

### Human IgG Specific, 550 Labeled

### 5ml Atomiser

### K8211

<b>Batch #</b>	<b>CDC0313A</b>
----------------	-----------------

<b>Full Description</b>	Goat anti-human IgG, polyclonal, F(ab)2, Dylight™ 550 conjugated. Sterile, azide free. In atomiser spray applicator
<b>Origin of Host Species</b>	Goat Serum
<b>Immunogen</b>	Human IgG
<b>Format</b>	F(ab)2 Conjugated to Dylight™ 550 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
<b>Specificity</b>	Human IgG
<b>Method of Purification</b>	Antigen affinity chromatography followed by pepsin digestion to obtain F(ab)2 fragments
<b>Appearance</b>	Clear pink solution
<b>Concentration</b>	2,000 U/ml
<b>Volume</b>	5.0 ml
<b>Total Amount</b>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>3.8 : 1</b>
<b>Buffer</b>	Phosphate Buffered Saline
<b>Preservatives/Stabilisers</b>	No preservatives
<b>Sterility</b>	0.2µm sterile filtered

Date of manufacture	20-March-2013
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**

*Christopher Lee*

Signed:

Date: 20 March 2013

Production Biologist