

# Certificate of Analysis

## CloneDetect

### Human IgG (Kappa Chain) Specific

### 650 Labeled

### K8251

<b>Batch #</b>	<b>CDS0714A</b>
<b>Full Description</b>	Goat anti-human IgG, (Kappa Chain), polyclonal, Dylight™ 650 conjugated. Sterile, azide free
<b>Origin of Host Species</b>	Goat Serum
<b>Immunogen</b>	Human IgG (Kappa Chain)
<b>Format</b>	IgG Conjugated to Dylight™ 650 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
<b>Specificity</b>	Human IgG (Kappa Chain)
<b>Method of Purification</b>	Affinity chromatography
<b>Appearance</b>	Clear blue solution
<b>Concentration</b>	10,000 U/ml
<b>Volume</b>	1.0 ml
<b>Total Amount</b>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>3.1</b>
<b>Buffer</b>	Phosphate Buffered Saline
<b>Preservatives/Stabilisers</b>	None
<b>Sterility</b>	0.2µm sterile filtered

<b>Date of manufacture</b>	<b>31-Jul-2014</b>
<b>Expiry Date</b>	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: 02-Sep-2014

*Christopher Lee*

**Production Biologist**