

# Certificate of Analysis

## CloneDetect

### Mouse IgG Specific, 549 Labeled

### K8230


<b>Batch #</b>	<b>CDG061A</b>
----------------	----------------

<b>Full Description</b>	Rabbit anti-mouse IgG, polyclonal, F(ab)2, Dylight™ 549 conjugated. Sterile, azide free
<b>Origin of Host Species</b>	Rabbit Serum
<b>Immunogen</b>	Mouse IgG
<b>Format</b>	F(ab)2 Conjugated to Dylight™ 549 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
<b>Specificity</b>	Mouse IgG (all sub-classes)
<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>	10,000 U/ml
<b>Volume</b>	1.0 ml
<b>Total Amount</b>	0.75 mg
<b>FITC : Protein Ratio</b>	<b>6.6 : 1.0</b>
<b>Buffer</b>	Phosphate Buffered Saline
<b>Preservatives/Stabilisers</b>	1% BSA (Sigma Product: A7906 Lot:038K0668)
<b>Sterility</b>	0.2µm sterile filtered

Date of manufacture	15/04/2009
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 -20°C**

Signed: 

Date: 09/07/10

**Production Supervisor**