

# Certificate of Analysis

## CloneDetect

### Human IgG (H+L) Specific, 650 Labelled

### K8212

<b>Batch #</b>	<b>CDW1212A</b>
----------------	-----------------

<b>Full Description</b>	Sheep anti-human IgG (H+L), polyclonal, DyLight™ 650 conjugated. Sterile, azide free
<b>Origin of Host Species</b>	USA Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision Inspected to confirm the absence of following: Foot and Mouth Disease; Rift Valley Fever; Rinderpest; Vesicular Stomatitis; Peste Des Petits Ruminants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie The sheep had no contact with animals of a lesser health.
<b>Immunogen</b>	Human IgG (H+L) purified from human serum
<b>Preparation</b>	Antiserum is prepared by immunising sheep with the above immunogen. Cross reactivity with bovine serum proteins has been minimised by insoluble adsorption. The antiserum is then conjugated with DyLight™ 650. Unreacted dye is removed by gel filtration.
<b>Format</b>	DyLight™ 650 conjugated IgG, liquid DyLight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
<b>Specificity</b>	The unconjugated immunoglobulin was shown to be specific by immunoelectrophoresis and double diffusion
<b>Method of Purification</b>	Affinity chromatography
<b>Functionality Test</b>	Tested and validated in functional assay on <i>ClonePix FL</i>
<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>3.3</b>
<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilisers</b>	None
<b>Sterility</b>	0.2µm sterile filtered
<b>Date of manufacture</b>	<b>12-December-2012</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: 26 December 2012

**Production Supervisor**