

Certificate of Analysis

CloneDetect

Human IgG (H+L) Specific, 650 Labelled K8212

Batch #	CDW0914A
----------------	-----------------

Full Description	Sheep anti-human IgG (H+L), polyclonal, DyLight™ 650 conjugated. Sterile, azide free
Origin of Host Species	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.
Immunogen	Human IgG (H+L) purified from human serum
Preparation	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.
Format	DyLight™ 650 conjugated IgG, liquid DyLight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
Specificity	The unconjugated immunoglobulin was shown to be specific by immunoelectrophoresis and double diffusion
Method of Purification	Affinity chromatography
Functionality Test	Tested and validated in functional assay on <i>ClonePix FL</i>
Appearance	Clear blue solution
Concentration	10,000 U/ml

Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	2.9
Buffer	Phosphate Buffered Saline pH7.2
Preservatives/Stabilisers	None
Sterility	0.2µm sterile filtered
Date of manufacture	09-Nov-2014
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: 20-Jan-2015

Christopher Lee

Production Biologist