

Certificate of Analysis

CloneDetect

Human IgG (Kappa Chain) Specific

550 Labeled

K8250

Batch #	CDQ0914A
Full Description	Sheep anti-human IgG, (Kappa Chain), polyclonal, Dylight™ 550 conjugated. Sterile, azide free, BSA free
Origin of Host Species	Sheep Serum
Immunogen	Human IgG (Kappa Chain)
Format	IgG Conjugated to Dylight™ 550 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
Specificity	Human IgG (Kappa Chain)
Method of Purification	Affinity chromatography
Appearance	Clear pink solution
Concentration	0.75 mg/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	4.0
Buffer	Phosphate Buffered Saline, pH 7.2
Preservatives/Stabilisers	None
Sterility	0.2µm sterile filtered

Date of manufacture	04-Sep-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 -20°C

Signed:

Date: 30-Sep-2014

Christopher Lee

Production Biologist