

Certificate of Analysis

CloneDetect

Human IgG (Kappa Chain) Specific

650 Labeled

K8251

Batch #	CDS0713A
Full Description	Goat anti-human IgG, (Kappa Chain), polyclonal, Dylight™ 650 conjugated. Sterile, azide free
Origin of Host Species	Goat Serum
Immunogen	Human IgG (Kappa Chain)
Format	IgG Conjugated to Dylight™ 650 – 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 blue solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	4.5
Buffer	Phosphate Buffered Saline
Preservatives/Stabilisers	None
Sterility	0.2µm sterile filtered

Date of manufacture	19 July 2013
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: 8 August 2013