

Certificate of Analysis

CloneDetect

Human IgG (Kappa Chain) Specific Fluorescein

K8252

Batch #	CDR1112A
Full Description	Sheep anti-Human kappa light chain (free and bound) Affinity Purified - Fluorescein, Azide free, BSA Free conjugate
Origin of Host Species	Sheep Serum
Immunogen	Human IgG (Kappa Chain)
Format	Fluorescein conjugated IgG – Liquid
Specificity	Human IgG (Kappa Chain)
Method of Purification	Affinity chromatography
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	3.7
Buffer	Phosphate Buffered Saline
Preservatives/Stabilisers	None
Sterility	0.2µm sterile filtered

Date of manufacture	08 November 2012
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: *Christopher Lee*

Date: 13 December 2012

Production Biologist