

Certificate of Analysis

CloneDetect

Human IgG (Kappa Chain) Specific Fluorescein

K8252

| | |
|----------------------------------|---|
| Batch # | CDR0515A |
| 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 | 4.5 |
| Buffer | Phosphate Buffered Saline |
| Preservatives/Stabilisers | None |
| Sterility | 0.2µm sterile filtered |

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 *Kieu Khiem Thi Nguyen*

Date: 17 December 2015

Production Biologist