

# Certificate of Analysis

## CloneDetect

### Human IgG (Kappa Chain) Specific Fluorescein

### K8252

<b>Batch #</b>	<b>CDR0717A</b>
<b>Full Description</b>	Sheep anti-Human kappa light chain (free and bound) Affinity Purified - Fluorescein, Azide free, BSA Free conjugate
<b>Origin of Host Species</b>	Sheep Serum
<b>Immunogen</b>	Human IgG (Kappa Chain)
<b>Format</b>	Fluorescein conjugated IgG – Liquid
<b>Specificity</b>	Human IgG (Kappa Chain)
<b>Method of Purification</b>	Affinity chromatography
<b>Appearance</b>	Clear yellow solution
<b>Concentration</b>	10,000 U/ml
<b>Volume</b>	1.0 ml
<b>Total Amount</b>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>5.3</b>
<b>Buffer</b>	Phosphate Buffered Saline
<b>Preservatives/Stabilisers</b>	None
<b>Sterility</b>	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: 21-MAR-2018

**Production Chemist III**