



# Certificate of Analysis

## CloneDetect Mouse IgG (H+L) Specific, Fluorescein 5ml Atomizer K8221

<b>Batch #</b>	<b>CDE0517A</b>
<b>Full Description</b>	Sheep anti-mouse IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile, azide free. In atomizer spray applicator
<b>Origin of Host Species</b>	USA Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision Inspected to confirm the absence of following: Foot and Mouth Disease; Rift Valley Fever; Rinderpest; Vesicular Stomatitis; Peste Des Petits Ruminants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie The sheep had no contact with animals of a lesser health.
<b>Immunogen</b>	Mouse IgG (H+L) purified from mouse serum
<b>Preparation</b>	Antiserum is prepared by immunising sheep with the above immunogen. Cross reactivity with bovine, human and rabbit serum proteins has been minimised by insoluble adsorption. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
<b>Format</b>	Fluorescein-conjugated IgG, liquid
<b>Specificity</b>	This product may give several arcs in the gamma region when tested by IEP against normal mouse serum (containing all four mouse IgG subclasses). Identity has been confirmed by double diffusion against mouse serum and anti-mouse immunoglobulin's of known specificity.
<b>Method of Purification</b>	Affinity chromatography
<b>Functionality Test</b>	This product contains K8220 diluted in PBS.
<b>Appearance</b>	Clear yellow solution
<b>Concentration</b>	2,000 U/ml
<b>Volume</b>	5.0 ml
<b>Total Amount</b>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>3.5</b>
<b>Buffer</b>	Phosphate Buffered Saline pH 7.2
<b>Preservatives/Stabilisers</b>	0.2% BSA (Sigma Product: <b>A3059</b> Lot <b>SLBK1938V</b> )
<b>Sterility</b>	0.2µm sterile filtered
<b>Expiry Date</b>	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 Nguyen*

Date: 23-May-2017

**Production Biologist**