

# Certificate of Analysis

## CloneDetect

### Human IgG (H+L) Specific, Fluorescein K8200

<b>Batch #</b>	<b>CDA0117A</b>
----------------	-----------------

<b>Full Description</b>	Sheep anti-human IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile, azide free
<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 Rumaninants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie The sheep had no contact with animals of a lesser health.
<b>Immunogen</b>	Human IgG (H+L) purified from human serum
<b>Preparation</b>	Antiserum is prepared by immunising sheep with the above immunogen. Cross reactivity with bovine 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>	The unconjugated immunoglobulin was shown to be specific by immunoelectrophoresis and double diffusion
<b>Method of Purification</b>	Affinity chromatography
<b>Viral Testing</b>	Negative. Product was found to be free of viral contaminants based on: qualified real-time PCR screening for rodent virus panels. Tests performed by Charles River (full report available on request).
<b>Mycoplasma Testing</b>	Negative. Product is assayed following a broth enrichment step using a <b>R&amp;D Testing for the Presence of Agar Cultivable and Noncultivable Mycoplasmas in Accordance with the United States Pharmacopoeia and the European Pharmacopoeia Guidelines GP-V611.19 assay</b> . Assay meets or exceeds the requirements of 21 CFR 610.30, 1993 PTC, and EP. Test performed by Charles River (report available on request).
<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>4.0</b>
<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilisers</b>	1% BSA (Manufacture <b>Boval</b> ; Catalog No: <b>LY-0082</b> ; Lot <b>LY-645</b> )
<b>Sterility</b>	0.2µm sterile filtered Product tested, no bacterial or fungal growth found after 72 hours incubation
<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: 10-MAR-2017

**Production Biologist**