



# Certificate of Analysis

## CloneDetect

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

### K8200

<b>Batch #</b>	CDA0818A
----------------	----------

<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.
<b>Immunogen</b>	Human IgG (H+L) purified from human serum
<b>Preparation</b>	Cross reactivity with bovine serum proteins has been minimised by affinity adsorption by binding to immobilised bovine proteins
<b>Format</b>	Affinity purified sheep immunoglobulin conjugated to fluorescein-5-EX, Succinimidyl Ester
<b>Specificity</b>	The unconjugated immunoglobulin was shown to be reactive to human serum and unreactive to mouse/rate/bovine serum by immunoelectrophoresis
<b>Method of Purification</b>	Affinity purification by absorption to immobilised human IgG
<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 Testing for Presence of Mycoplasma Using Broth Culture Growth Enrichment and RT-PCR Detection assay GP.V900.01. 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>3.5</b>
<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilizers</b>	1% BSA (Manufacture: Proliant; Catalog #: 68100 ; Lot # <b>BB71160101</b> )
<b>Sterility</b>	0.2µm sterile filtered.
<b>Expiry Date</b>	12 months from receipt of product
<b>Storage</b>	2 – 8°C

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

**For research use only**

Signed: *Kieu-Khiem Nguyen*

Date: 03-SEP-2019

**Production Biologist**