

# Certificate of Analysis

## CloneDetect

### Human IgG (Fc) Specific, Fluorescein K8205

<b>Batch #</b>	<b>CDB1212A</b>
----------------	-----------------

<b>Full Description</b>	Sheep anti-human IgG (Fc), 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 Ruminants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie The sheep had no contact with animals of a lesser health.
<b>Immunogen</b>	Human IgG (Fc) purified from human serum and digested with papain. Fc fraction purified and shown to be homogenous by SDS-PAGE
<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>	Polyclonal sheep immunoglobulin, fluorescein conjugated, liquid
<b>Specificity</b>	Cross reactivity with bovine, mouse and rat serum proteins minimised by insoluble adsorption. Shown to be specific by immunoelectrophoresis and double diffusion
<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>2.8</b>

<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilisers</b>	1% BSA (Sigma Product: A3059 Lot: 120M1885V)
<b>Sterility</b>	0.2µm sterile filtered Product tested, no bacterial or fungal growth found after 72 hours incubation
<b>Date of manufacture</b>	<b>12-Dec-2012</b>
<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°C – 8°C**

Signed:

Date: 20-Dec-2012

*Christopher Lee*

**Production Chemist**