

Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific, Fluorescein

5ml Atomiser

K8206

Batch #	CDB1212A
----------------	-----------------

Full Description	Sheep anti-human IgG (Fc), polyclonal, Fluorescein-conjugated. Sterile, azide free. In atomiser spray applicator
Origin of Host Species	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.
Immunogen	Human IgG (Fc) purified from human serum and digested with papain. Fc fraction purified and shown to be homogenous by SDS-PAGE
Preparation	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.
Format	Polyclonal sheep immunoglobulin, fluorescein conjugated, liquid
Specificity	Cross reactivity with bovine, mouse and rat serum proteins minimised by insoluble adsorption. Shown to be specific by immunoelectrophoresis and double diffusion
Method of Purification	Affinity chromatography
Appearance	Clear yellow solution
Concentration	2,000 U/ml
Volume	5.0 ml
Total Amount	0.75 mg

Dye : Protein Ratio	2.8
Buffer	Phosphate Buffered Saline pH7.2
Preservatives/Stabilisers	0.2% BSA (Sigma Product: A3059 Lot: 120M1885V)
Sterility	0.2µm sterile filtered Product tested, no bacterial or fungal growth found after 72 hours incubation
Date of manufacture	12-Dec-2012
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°C – 8°C

Signed:

Date: 20-Dec-2012

Christopher Lee

Production Chemist