



Certificate of Analysis

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

Batch #	CDE0315A
Full Description	Sheep anti-mouse IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile, azide free. In atomizer 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	Mouse IgG (H+L) purified from mouse serum
Preparation	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.
Format	Fluorescein-conjugated IgG, liquid
Specificity	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.
Method of Purification	Affinity chromatography
Functionality Test	This product contains K8220 diluted in PBS.
Appearance	Clear yellow solution
Concentration	2,000 U/ml
Volume	5.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	2.9
Buffer	Phosphate Buffered Saline pH 7.2
Preservatives/Stabilisers	0.2% BSA (Sigma Product: A3059 Lot SLBD5581V)
Sterility	0.2µm sterile filtered
Date of manufacture	11-Mar-2015
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 – 8°C

Signed: Christopher Lee
Production Biologist

Date: 11-Mar-2015