

Certificate of Analysis

CloneDetect

Mouse IgG (Fc) Specific, Fluorescein K8225

Batch #	CDF0914A
----------------	-----------------

Full Description	Sheep anti-mouse IgG (Fc), polyclonal, Fluorescein-conjugated. Sterile, azide free
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 (Fc). IgG (all subclasses) purified from mouse serum and digested with papain. Fc fraction purified and shown to be >95% pure by SDS-PAGE
Preparation	Antiserum is prepared by immunising sheep with the above immunogen. Immunoglobulin affinity purified and conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
Format	Polyclonal sheep immunoglobulin, fluorescein-conjugated, liquid
Specificity	Cross reactivity with bovine, human and rabbit serum proteins minimised by insoluble adsorption. Shows <5% cross-reactivity with IgA and IgM by ELISA. Identity confirmed by double diffusion against mouse serum and anti-mouse IgG of known specificity. Gives several arcs when tested by IEP against normal mouse serum (mouse IgG subclasses).
Method of Purification	Affinity chromatography
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg

Dye : Protein Ratio	2.8
Buffer	Phosphate Buffered Saline pH 7.2
Preservatives/Stabilisers	1% BSA (Sigma Product: A3059 Lot SLBD5581V)
Sterility	0.2µm sterile filtered
Date of manufacture	08-SEP-2014
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: 

Date: 04-Nov-2014

Production Biologist