



Certificate of Analysis

CloneDetect

Mouse IgG (Fc) Specific, Fluorescein

K8225

| | |
|----------------|-----------------|
| Batch # | CDF0117A |
|----------------|-----------------|

| | |
|-------------------------------|--|
| 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 | 3.2 |
| Buffer | Phosphate Buffered Saline pH 7.2 |
| Preservatives/Stabilisers | 1% BSA (Product Code: LY-0082 ; Lot LY-645) |
| Sterility | 0.2µm sterile filtered |
| 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:

Kieu Khiem Nguyen

Date: 17- FEB-2017

Production Biologist