



# Certificate of Analysis

## CloneDetect

### Mouse IgG (Fc) Specific, Fluorescein

## K8225

|                |                 |
|----------------|-----------------|
| <b>Batch #</b> | <b>CDF0119B</b> |
|----------------|-----------------|

|                               |  |
|-------------------------------|--|
| <b>Full Description</b>       | Sheep anti-mouse IgG (Fc), polyclonal, Fluorescein-conjugated. Sterile, azide free   |
| <b>Origin of Host Species</b> | USA<br>Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision<br>Inspected to confirm the absence of following:<br>Foot and Mouth Disease; Rift Valley Fever; Rinderpest; Vesicular Stomatitis; Peste Des Petits Ruminants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie<br>The sheep had no contact with animals of a lesser health. |
| <b>Immunogen</b>              | 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  |
| <b>Preparation</b>            | Antiserum is prepared by immunising sheep with the above immunogen. Immunoglobulin affinity purified and 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, 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).  |
| <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>4.9</b>  |
| <b>Buffer</b>                    | Phosphate Buffered Saline pH 7.2                          |
| <b>Preservatives/Stabilisers</b> | 1% BSA ( <b>Boval Product Code: LY-0082; Lot LY-749</b> ) |
| <b>Sterility</b>                 | 0.2µm sterile filtered                                    |
| <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 – 8°C**

Signed:

*Kieu Khiem Nguyen*

Date: 26-JUN-2019

**Reagent Biologist**