



Certificate of Analysis

CloneDetect

Mouse IgG (Fc)_γ Specific, Fluorescein K8425

| | |
|----------------|---------------|
| Batch # | 170218 |
|----------------|---------------|

| | |
|-------------------------------|---|
| Full Description | Goat anti-mouse IgG, Fc _γ fragment specific, polyclonal, Fluorescein (FITC)-conjugated. Sterile, azide free |
| Origin of Host Species | USA Goat kept at a USDA registered technical plant (registration # 23-R-0114) under regular veterinary supervision and inspection. Raw materials are collected from animals veterinary-certified clinically free of signs of bovine spongiform encephalopathy (BSE), transmissible spongiform encephalopathy (TSE), foot and mouth disease (FMD). The goat had no contact with animals of a lesser health and have remained healthy during raw materials collection. |
| Immunogen | Mouse IgG (Fc) |
| Preparation | Antiserum is prepared by immunizing goat with the above immunogen. Immunoglobulin affinity purified and conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration. |
| Format | Polyclonal goat immunoglobulin, fluorescein-conjugated, liquid |
| Specificity | Based on Immuno-electrophoresis and/or ELISA, the product reacts with the Fc portion of mouse IgG heavy chain but not with the Fab portion of mouse immunoglobulins. No cross-reactivity was detected against mouse IgM or against non-immunoglobulin serum proteins. |
| 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 | 3.5 |
| Buffer | Phosphate Buffered Saline pH 7.6 |
| Preservatives/Stabilizers | 1.5% Bovine Serum Albumin (IgG-Free, Protease-Free) |
| Sterility | 0.2µm sterile filtered |
| Expiry Date | 6 months from shipment of product |

This product has been manufactured to ISO 9001:2015 Quality Assurance guidelines

For research use only Store at 2 – 8°C, avoid direct light

Signed: *Kieu Khiem Nguyen*

Date: 19-MAR-2024

Production Chemist