



Certificate of Analysis

CloneDetect

Mouse IgG (Fc)_γ Specific, Fluorescein K8425

Batch #	158607
----------------	---------------

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

Document Revised on September 2, 2022 to correct "Expiry Date" from "9 months from shipment of product" to "6 months from shipment of product".

Dye : Protein Ratio	3.9
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: *Erwin Ramirez*

Date: 24-Jan-2022

Production Chemist

Document Revised on September 2, 2022 to correct "Expiry Date" from "9 months from shipment of product" to "6 months from shipment of product".