



Certificate of Analysis

CloneDetect

Mouse IgG (H+L) Specific, Fluorescein K8420

Batch #	165238
----------------	---------------

Full Description	Goat anti-mouse IgG (H+L), 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 purified from mouse serum
Preparation	Antiserum is prepared by immunizing goat with the above immunogen. Cross reactivity with bovine, human, horse, swine and rabbit serum proteins has been minimized by insoluble adsorption. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
Format	Polyclonal goat immunoglobulin, fluorescein-conjugated, liquid
Specificity	This product reacts with whole molecule mouse IgG. It may give several arcs in the gamma region when tested by IEP against normal mouse serum. Identity has been confirmed by immuno-electrophoresis and/or ELISA against mouse serum and anti-mouse immunoglobulins with known specificity.
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	3.4
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: 18-MAY-2023

Production Chemist