



# Certificate of Analysis

## CloneDetect

### Mouse IgG (Fc)<sub>γ</sub> Specific, Fluorescein K8425

<b>Batch #</b>	<b>161223</b>
----------------	---------------

<b>Full Description</b>	Goat anti-mouse IgG, Fc <sub>γ</sub> fragment specific, polyclonal, Fluorescein (FITC)-conjugated. Sterile, azide free
<b>Origin of Host Species</b>	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.
<b>Immunogen</b>	Mouse IgG (Fc)
<b>Preparation</b>	Antiserum is prepared by immunizing goat with the above immunogen. Immunoglobulin affinity purified and conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
<b>Format</b>	Polyclonal goat immunoglobulin, fluorescein-conjugated, liquid
<b>Specificity</b>	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.
<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

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

<b>Dye : Protein Ratio</b>	<b>3.2</b>
<b>Buffer</b>	Phosphate Buffered Saline pH 7.6
<b>Preservatives/Stabilizers</b>	1.5% Bovine Serum Albumin (IgG-Free, Protease-Free)
<b>Sterility</b>	0.2µm sterile filtered
<b>Expiry Date</b>	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: 03-AUG-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".