



# Certificate of Analysis

## CloneDetect

### Mouse IgG (H+L) Specific, Fluorescein K8420

<b>Batch #</b>	<b>158188</b>
----------------	---------------

<b>Full Description</b>	Goat anti-mouse IgG (H+L), 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 purified from mouse serum
<b>Preparation</b>	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.
<b>Format</b>	Polyclonal goat immunoglobulin, fluorescein-conjugated, liquid
<b>Specificity</b>	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.
<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>3.9</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: Ervin Ramirez

Date: 06-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".