

# Certificate of Analysis

## CloneDetect

### Human IgG (H+L) Specific, Fluorescein K8400

<b>Batch #</b>	<b>159353</b>
----------------	---------------

<b>Full Description</b>	Goat anti-human IgG (H+L), polyclonal, FITC Fluorescein-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>	Human IgG (H+L) purified from human serum
<b>Preparation</b>	Antiserum is prepared by immunizing goat with the above immunogen. Cross reactivity with bovine, horse, mouse serum proteins has been minimized by insoluble adsorption. The purified immunoglobulin is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
<b>Format</b>	Affinity purified goat immunoglobulin conjugated to fluorescein-5-isothiocyanate
<b>Specificity</b>	This product reacts with whole molecule human IgG. It may also react with the light chains of other human immunoglobins. No antibody was detected against non-immunoglobulin serum proteins by immunoelectrophoresis and/or ELISA.
<b>Method of Purification</b>	Immunoaffinity chromatography using antigens coupled to agarose beads
<b>Purified Antibody/Starting Materials Identification ID</b>	<b>159353</b>
<b>Viral Testing</b>	Negative. Product was found to be free of viral contaminants based on: qualified real-time PCR screening for comprehensive rodent virus panels. Tests performed by Charles River (report available on request).

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>Mycoplasma Testing</b>	Negative. Product is assayed following a broth enrichment step using a Testing for Presence of Mycoplasma Using Broth Culture Growth Enrichment and RT-PCR Detection assay GP.V900.01. Assay meets or exceeds the requirements of 21 CFR 610.30, 1993 PTC, and EP. Test performed by Charles River (report available on request).
<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
<b>Dye : Protein Ratio</b>	➤ <b>3.5</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
<b>Storage</b>	2 – 8°C

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

**For research use only**

Signed: *Kieu-Khiem Nguyen*

Date: 23-AUG-2022

**Production Biologist**