



# Certificate of Analysis

## CloneDetect

### Human IgG (Fc) Specific

### Recombinant, Fluorescein

## K8295

<b>Batch #</b>	<b>CDU0417A</b>
<b>Full Description</b>	Mouse anti-human IgG (Fc), recombinant IgG1. Fluorescein-conjugated, sterile, azide free. Does not contain products of animal origin (for further information refer to Certificate of Origin).
<b>Origin of Host Species</b>	Chinese Hamster Ovary (CHO) cells, grown in animal-free conditions.
<b>Format</b>	Fluorescein conjugated (NHS Ester) IgG, liquid
<b>Specificity</b>	Human IgG Fc (all subclasses)
<b>Purified Mouse IgG1 Antibody</b>	MCA647-MK1A6 Batch 16-957
<b>Method of Purification</b>	Protein A 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
<b>Dye : Protein Ratio</b>	4.3
<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilisers</b>	None
<b>Viral Testing</b>	Negative. Bulk harvest assessed for presence of viral contamination using a mouse/rat comprehensive PCR panel. Test performed by Charles River (further details and full report available on request).
<b>Mycoplasma Testing</b>	Negative. Bulk harvest tested by RT-PCR. Test performed by Charles River (further details and full report available on request).
<b>Endotoxin Testing</b>	6.49 EU/ml by LAL method Test performed by BioReliance (further details and full report available on request).
<b>Sterility</b>	0.2µm sterile filtered
<b>Expiry Date</b>	12 months from receipt of product

This product has been manufactured to ISO9001:2015 Quality Assurance guidelines

**For research use only    Store at 2 – 8°C**

Signed: *Kieu-Khiem Nguyen*

Date: 21 JUN-2017



**Production Biologist**

# Certificate of Origin

## CloneDetect

### Human IgG (Fc) Specific

### Recombinant, Fluorescein

## K8295

K8295 is a recombinant fluorescein-conjugated monoclonal mouse IgG1 against human IgG produced using CHO (Chinese Hamster Ovary) cells. The final antibody is formulated in sterile PBS free of azide. To the best of our knowledge and according to information obtained from our suppliers, K8295 contains no raw materials of animal origin, nor are any materials of animal origin used in the manufacture of the antibody, and the product has not been exposed to BSE/ TSE (bovine/ transmissible spongiform encephalopathies).

Signed: Signed: *Kieu-Khiem Nguyen*

Name: Kieu-Khiem Thi Nguyen

Position: Production Biologist

Date: 21-JUN-2017