



Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific

Recombinant, Fluorescein

K8295

Batch #	CDU0618A
Full Description	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).
Origin of Host Species	Chinese Hamster Ovary (CHO) cells, grown in animal-free conditions.
Format	Fluorescein conjugated (NHS Ester) IgG, liquid
Specificity	Human IgG Fc (all subclasses)
Purified Mouse IgG1 Antibody	MCA647-MK1A6 Batch 18-1057
Method of Purification	Protein A affinity chromatography
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	4.0
Buffer	Phosphate Buffered Saline pH7.2
Preservatives/Stabilisers	None
Viral Testing	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).
Mycoplasma Testing	Negative. Bulk harvest tested by RT-PCR. Test performed by Charles River (further details and full report available on request).
Endotoxin Testing	< 0.5 EU/ml by LAL method Test performed by BioReliance (further details and full report available on request).
Sterility	0.2µm sterile filtered
Expiry Date	12 months from receipt of product

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

For research use only Store at 2 – 8°C

Signed: *Kieu Khiem Nguyen*

Date: 31- JUL- 2018

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: *Kieu Khiem Nguyen*

Name: Kieu Khiem Nguyen

Position: Chemist III

Date: 31- JUL- 2018