



Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific Recombinant, Fluorescein

K8295

| | |
|-------------------------------------|--|
| Batch # | CDU1118A |
| 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 18-1121 |
| 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 | ➤ 2.8 / 4.0 |
| Buffer | 0.145M PBS pH 7.2 |
| Preservatives/Stabilisers | No BSA; No Azide |
| 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 | 15.641 EU/ml by Kinetic-Chromogenic LAL method. Test performed by Charles River Laboratory (further details and full report available on request). |



| | |
|--------------------|-----------------------------------|
| Sterility | 0.2µm sterile filtered |
| Expiry Date | 12 months from receipt of product |
| Storage | 2 – 8°C |

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

For research use only

Signed: *Kieu Khiem Nguyen*

Date: 25 - OCT - 2019

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: 25- OCT -2019