



Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific Recombinant, Fluorescein

K8495

Batch #	1033109
---------	---------

Full Description	Mouse anti-human IgG (Fc), recombinant IgG1. FITC Fluorescein-conjugated, monoclonal, 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 origin-free conditions
Format	Fluorescein conjugated (NHS Ester) IgG, liquid
Specificity	Human IgG Fc (all subclasses)
Purified Mouse IgG1 Antibody	MCA647-MK1A6 Batch 19-0422
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
Buffer	0.145M PBS pH 7.2
Preservatives/Stabilizers	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	<0.05 EU/ml by LAL method Test performed by CRL (further details and full report available on request).
Sterility	0.2µm sterile filtered



Expiry Date	6 months from shipment of product
Storage	2 – 8°C

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

For research use only

Signed: *Kieu Khiem nguyen*

Date: 28-NOV- 2023

Production Biologist



Certificate of Origin

CloneDetect

Human IgG (Fc) Specific Recombinant, Fluorescein K8495

K8495 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, K8495 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

Date: 28-NOV 2023