



Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific, Fluorescein

K8205

Batch #	CDB0818A
----------------	-----------------

Full Description	Sheep anti-human IgG (Fc), polyclonal, Fluorescein-conjugated. Sterile, azide free
Origin of Host Species	USA Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision.
Immunogen	Human IgG (Fc) purified from human serum and digested with papain. Fc fraction purified and shown to be homogenous by SDS-PAGE
Preparation	Antiserum is prepared by immunising sheep with the above immunogen. Cross reactivity with bovine serum proteins has been minimised by insoluble adsorption. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
Format	Polyclonal sheep immunoglobulin, fluorescein conjugated, liquid
Specificity	The unconjugated immunoglobulin was shown to be reactive to human serum and unreactive to mouse/rate/bovine serum by immune-electrophoresis
Method of Purification	Affinity purification by absorption to immobilised human IgG
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	> 2.8 / 4.3
Buffer	Phosphate Buffered Saline pH7.2
Preservatives/Stabilisers	1% BSA (Proliant Cat # 68100; Lot: # BB71160101)
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: 12 – OCT- 2018

Production Biologist