



Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific, Fluorescein

K8205

| | |
|----------------|-----------------|
| Batch # | CDB0719A |
|----------------|-----------------|

| | |
|----------------------------------|---|
| 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 / 3.8 |
| 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: 15 – APR- 2020

Production Biologist