

Certificate of Analysis

CloneDetect

Human IgG Specific, 649 Labeled

5ml Atomiser

K8213

Batch #	CDW0516A
----------------	-----------------

Full Description	Goat anti-human IgG, polyclonal, F(ab)2, Dylight™ 649 conjugated. Sterile, azide free. In atomiser spray applicator
Origin of Host Species	Goat Serum
Immunogen	Human IgG
Format	F(ab)2 Conjugated to Dylight™ 649 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
Specificity	Human IgG
Method of Purification	Antigen affinity chromatography followed by pepsin digestion to obtain F(ab)2 fragments
Appearance	Clear blue solution
Concentration	2,000 U/ml
Volume	5.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	3.3
Buffer	Phosphate Buffered Saline
Sterility	0.2µm sterile filtered

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

For research use only Store at -20°C

Signed: *Kieu-Khiem Nguyen*

Date: 27 MAY 2016

Production Chemist