

Certificate of Analysis

CloneDetect

Mouse IgG Specific, 650 Labeled

K8233

Batch #	CDX0516A
----------------	-----------------

Full Description	Rabbit anti-mouse IgG, polyclonal, F(ab)2, Dylight™ 650 conjugated. Sterile, azide free
Origin of Host Species	Rabbit Serum
Immunogen	Mouse IgG
Format	F(ab)2 Conjugated to Dylight™ 650 – Liquid Dylight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries
Specificity	Mouse IgG (all sub-classes)
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
FITC : Protein Ratio	4.0 : 1.0
Buffer	Phosphate Buffered Saline
Preservatives/Stabilisers	1% BSA (Sigma Product:A7906 Lot:069K1653)
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 Supervisor