
Certificate of Analysis

K8255
CloneDetect
Anti-Human
Detection Agent
IgM Specific
FITC Labelled

Batch #	CDP0509A
Full Description	Goat anti Human IgM FITC: Azide free
Origin of Host Species	Goat Serum
Immunogen	Human IgM
Purification method	Affinity chromatography
Format	IgG FITC conjugated – liquid
Specificity	Human
Appearance	Clear yellow solution
Concentration	10,000 U/mL
Volume	1.0 mL
Total Amount	0.75mg
FITC : Protein Ratio	9.2 : 1.0
Buffer	Phosphate Buffered Saline
Stabilisers	1% BSA (Sigma Product:A7906 Lot:038K0668)
Date of manufacture	April 2009
Expiry Date	12 months from receipt of product

This product has been manufactured to ISO9001 Quality Assurance guidelines

For research use only Store at -20°C

Signed  Date 11/05/09

Reagents Laboratory Team Leader