
Certificate of Analysis

K8253
CloneDetect Anti-Human
Detection Agent
IgM Specific
549 Labelled

Genetix

BATCH #	CDJ1208/84805
Full Description	Goat anti human IgM : Dylight™ 549 labelled
Origin of Host Species	Goat Serum
Immunogen	Human IgM
Format	IgG Conjugated to Dylight 549 DyLight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries.
Specificity	Human
Method of Purification	Affinity Chromatography
Appearance	Clear, pink liquid
Concentration	10 000 U/mL
Volume	1 mL
Total Amount	0.75 mg
549 : Protein Ratio	4.3 : 1.0
Buffer	Phosphate Buffered Saline
Stabilisers/ Preservatives	1% BSA
Sterility	0.2µm Sterile Filtered
Date of manufacture	18/11/08

This product has been manufactured to ISO9001:2000
For research use only Store at -20°C

Signed  **Date** 3/12/08

Reagents Laboratory Team Leader