
Certificate of Analysis

K8254 CloneDetect Anti-
Human Detection Agent
IgM Specific 649 Labelled

Batch # **CDL0208/72291**

Full Description Goat anti human igM : DyLight™¹ 649 labelled

Origin of Host Species Goat Serum

Immunogen Human IgM

Preparation Antisera to human IgM were raised by repeated immunisation of goats with highly purified antigen. Purified igG was prepared from whole serum by affinity chromatography.

Format DyLight™¹649 conjugated IgG

Specificity Human IgM

Method of Purification Affinity chromatography

Over /

Appearance	Clear blue solution solution
Concentration	10,000 U/mL
Volume	1.0 mL
Total Amount	0.75 mg
649 : Protein Ratio	5.3:1.0
Buffer	Phosphate Buffered Saline
Stabilisers	1% BSA (certificate available on request)
Preservatives	None
Sterility	0.2µm sterile filtered
Date of manufacture	21/01/08

This product has been manufactured to ISO9000 & ISO13485 Quality Assurance guidelines.

For research use only. Store at -20°C

Signed  _____ Date _____

Reagents Laboratory Manager

¹DyLight™ is a trademark of Thermo Fisher Scientific Inc. and its subsidiaries.

Europe

Genetix
Queensway, New Milton,
Hampshire, BH25 5AN, UK
Tel +44 (0) 1425 624758
Fax +44 (0) 1425 624747

Genetix

Humboldtstr. 12,
D-85609 Gernach, Munich, Germany
Tel +49 (0) 89 844 80 275/276
Fax +49 (0) 89 844 80 110

USA

Genetix
56 Roland Street,
Boston, MA 02129, USA
Tel +1 (617) 776 6100
Fax +1 (248) 522 7489

Genetix

120 Baytech Drive,
San José, CA 95134-2302, USA
Tel +1 (408) 719 6400
Fax +1 (408) 719 6401