

Certificate of Analysis

CloneDetect

Rat IgG (H+L) Specific, Fluorescein K8240

Batch #	CDI1115A
----------------	-----------------

Full Description	Sheep anti-rat IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile
Origin of Host Species	<p>USA</p> <p>Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision</p> <p>Inspected to confirm the absence of following:</p> <p>Foot and Mouth Disease; Rift Valley Fever; Rinderpest; Vesicular Stomatitis; Peste Des Petits Rumaninants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie</p> <p>The sheep had no contact with animals of a lesser health.</p>
Immunogen	Rat IgG (H+L) purified from rat serum
Preparation	Antiserum is prepared by immunising sheep with the above immunogen. Cross reactivity with bovine, human and rabbit serum proteins has been minimised by insoluble adsorption. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
Format	FITC-conjugated IgG, liquid
Specificity	This product may give several arcs in the gamma region when tested by IEP against rat serum (rat IgG subtypes and IgM). Identity has been confirmed by double diffusion against rat serum and anti-rat immunoglobulin's of known specificity.
Method of Purification	Affinity chromatography
Appearance	Clear yellow solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	1.0 mg

Dye : Protein Ratio	4.1
Buffer	Phosphate Buffered Saline pH 7.2
Preservatives/Stabilisers	1% BSA (Sigma Product: A3059 Lot SLBM7676V)
Sterility	0.2µm sterile filtered
Expiry Date	12 months from receipt of product

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

For research use only Store at 2 – 8°C

Signed:

Date: 25-FEB-2016

Kieu-Khiem Nguyen

Production Biologist