

# Certificate of Analysis

## CloneDetect

### Rat IgG (H+L) Specific, Fluorescein K8240

<b>Batch #</b>	<b>CDI0120A</b>
----------------	-----------------

<b>Full Description</b>	Sheep anti-rat IgG (H+L), polyclonal, Fluorescein-conjugated. Sterile
<b>Origin of Host Species</b>	<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 Ruminants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie</p> <p>The sheep had no contact with animals of a lesser health.</p>
<b>Immunogen</b>	Rat IgG (H+L) purified from rat serum
<b>Preparation</b>	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.
<b>Format</b>	FITC-conjugated IgG, liquid
<b>Specificity</b>	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.
<b>Method of Purification</b>	Affinity chromatography
<b>Appearance</b>	Clear yellow solution
<b>Concentration</b>	10,000 U/ml
<b>Volume</b>	1.0 ml
<b>Total Amount</b>	1.0 mg

<b>Dye : Protein Ratio</b>	<b>3.2</b>
<b>Buffer</b>	Phosphate Buffered Saline pH 7.2
<b>Preservatives/Stabilisers</b>	1% BSA ( <b>Proliant</b> Catalog #: 68100; lot # BB71870101)
<b>Sterility</b>	0.2µm sterile filtered
<b>Expiry Date</b>	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: 15-APR-2020

*Kieu-Khiem Nguyen*

**Production Biologist**