



# Certificate of Analysis

## CloneDetect

### Sheep IgG (H+L) Specific, Fluorescein

## K8246

<b>Batch #</b>	<b>CDK0418A</b>
<b>Full Description</b>	Donkey anti-sheep IgG (H+L), polyclonal, fluorescein-conjugated. Sterile, azide-free
<b>Origin of Host Species</b>	Donkey
<b>Immunogen</b>	Sheep IgG (H+L) purified from sheep serum
<b>Preparation</b>	Antiserum to sheep IgG was raised by repeated immunisation of donkeys with highly purified antigen. Purified IgG was prepared by affinity chromatography. The antiserum is then conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
<b>Format</b>	Donkey IgG (H+L) Fluorescein conjugated IgG, liquid
<b>Specificity</b>	The unconjugated immunoglobulin was shown to be specific by immunoelectrophoresis and double diffusion.
<b>Method of Purification</b>	Affinity chromatography
<b>Appearance</b>	Clear colourless solution
<b>Concentration</b>	10,000 U/ml
<b>Volume</b>	1.0 ml
<b>Total Amount</b>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>4.0</b>
<b>Buffer</b>	Phosphate Buffered Saline pH 7.2
<b>Sterility</b>	0.2µm sterile filtered Product tested, no bacterial or fungal growth found after 72 hours incubation
<b>Preservatives/Stabilisers</b>	1% BSA (Proliant Product: Cat #: 68100; Lot: BB70450101)
<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: *Kieu-Khiem Nguyen*

Date: 09-MAY – 2018

Production Biologist