



Certificate of Analysis

CloneDetect

Sheep IgG (H+L) Specific, Fluorescein

K8246

Batch #	CDK0419A
Full Description	Donkey anti-sheep IgG (H+L), polyclonal, fluorescein-conjugated. Sterile, azide-free
Origin of Host Species	Donkey
Immunogen	Sheep IgG (H+L) purified from sheep serum
Preparation	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.
Format	Donkey IgG (H+L) Fluorescein conjugated IgG, liquid
Specificity	The unconjugated immunoglobulin was shown to be specific by immunoelectrophoresis and double diffusion.
Method of Purification	Affinity chromatography
Appearance	Clear colourless solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Dye : Protein Ratio	3.5
Buffer	Phosphate Buffered Saline pH 7.2
Sterility	0.2µm sterile filtered Product tested, no bacterial or fungal growth found after 72 hours incubation
Preservatives/Stabilisers	1% BSA (Proliant Product: Cat #: 68100; Lot: BB71870101)
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: *Kieu-Khiem Nguyen*

Date: 15- MAY- 2019

Production Biologist