

# Certificate of Analysis

## CloneDetect

### Mouse IgG (Fc) Specific, Fluorescein 5ml Atomiser K8226

<b>Batch #</b>	<b>CDF1219A</b>
----------------	-----------------

<b>Full Description</b>	Sheep anti-mouse IgG (Fc), polyclonal, Fluorescein-conjugated. Sterile, azide free. In atomiser spray applicator
<b>Origin of Host Species</b>	USA Sheep kept at a registered technical plant (1774/2002 registration # VT-TEC-0002) under regular veterinary supervision Inspected to confirm the absence of following: Foot and Mouth Disease; Rift Valley Fever; Rinderpest; Vesicular Stomatitis; Peste Des Petits Rumaninants; Blue Tongue; Epizootic Hemorrhagic Disease; Scrapie The sheep had no contact with animals of a lesser health.
<b>Immunogen</b>	Mouse IgG (Fc). IgG (all subclasses) purified from mouse serum and digested with papain. Fc fraction purified and shown to be >95% pure by SDS-PAGE
<b>Preparation</b>	Antiserum is prepared by immunising sheep with the above immunogen. Immunoglobulin affinity purified and conjugated with fluorescein. Unreacted fluorochrome is removed by gel filtration.
<b>Format</b>	Polyclonal sheep immunoglobulin, fluorescein-conjugated, liquid
<b>Specificity</b>	Cross reactivity with bovine, human and rabbit serum proteins minimised by insoluble adsorption. Shows <5% cross-reactivity with IgA and IgM by ELISA. Identity confirmed by double diffusion against mouse serum and anti-mouse IgG of known specificity. Gives several arcs when tested by IEP against normal mouse serum (mouse IgG subclasses).
<b>Method of Purification</b>	Affinity chromatography
<b>Functionality Test</b>	This product contains K8225 diluted in PBS, the K8225 component is tested and validated in functional assay on <i>ClonePix FL</i>
<b>Appearance</b>	Clear yellow solution
<b>Concentration</b>	2,000 U/ml
<b>Volume</b>	5.0 ml
<b>Total Amount</b>	0.75 mg

<b>Dye : Protein Ratio</b>	<b>3.3</b>
<b>Buffer</b>	Phosphate Buffered Saline pH 7.2
<b>Preservatives/Stabilisers</b>	0.2% BSA ( <b>Proliant Biologicals Cat #: 68100; Lot BB71870101</b> )
<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:** *Kieu-Khiem Nguyen*

**Date:** 13-MAR-2020

**Production Biologist**