

# Certificate of Analysis

## CloneDetect

### Human IgG (Fc) Specific Recombinant, Fluorescein K8295

<b>Batch #</b>	<b>CDU0513A</b>
----------------	-----------------

<b>Full Description</b>	Mouse anti-human IgG (Fc), recombinant IgG1. Fluorescein-conjugated, sterile, azide free. Does not contain products of animal origin (for further information refer to Certificate of Origin).
<b>Origin of Host Species</b>	Chinese Hamster Ovary (CHO) cells, grown in animal-free conditions.
<b>Format</b>	Fluorescein conjugated (NHS Ester) IgG, liquid
<b>Specificity</b>	Human IgG Fc (all subclasses)
<b>Method of Purification</b>	Protein A 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>	0.75 mg
<b>Dye : Protein Ratio</b>	<b>3.3</b>
<b>Buffer</b>	Phosphate Buffered Saline pH7.2
<b>Preservatives/Stabilisers</b>	None
<b>Viral Testing</b>	Negative. Bulk harvest assessed for presence of viral contamination using a mouse/rat comprehensive PCR panel. Test performed by Charles River (further details and full report available on request).
<b>Mycoplasma Testing</b>	Negative. Bulk harvest tested by RT-PCR. Test performed by Charles River (further details and full report available on request).
<b>Endotoxin Testing</b>	< <b>8.9</b> Eu/ml by LAL method Test performed by BioReliance (further details and full report available on request).
<b>Sterility</b>	0.2µm sterile filtered

Date of manufacture	09 May 2013
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: 12 June 2013

**Production Supervisor**

# Certificate of Origin

**CloneDetect**

**Human IgG (Fc) Specific  
Recombinant, Fluorescein  
K8295**

K8295 is a recombinant fluorescein-conjugated monoclonal mouse IgG1 against human IgG produced using CHO (Chinese Hamster Ovary) cells. The final antibody is formulated in sterile PBS free of azide. To the best of our knowledge and according to information obtained from our suppliers, K8295 contains no raw materials of animal origin, nor are any materials of animal origin used in the manufacture of the antibody, and the product has not been exposed to BSE/ TSE (bovine/ transmissible spongiform encephalopathies).



Signed:

Name: Meguel Varda

Position: Production Supervisor

Date: 4 June 2013