

Certificate of Analysis

CloneDetect

Human IgG (Fc) Specific CIF (Complex Initiation Factor)

K8215

Batch #	CDD0413A
----------------	-----------------

Full Description	Sheep anti-human IgG (H+L), polyclonal, unconjugated. Sterile, azide free
Origin of Host Species	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.
Immunogen	Human IgG (H+L) purified from human serum
Format	Non-conjugated IgG, liquid
Specificity	This product gives several arcs in the gamma region by IEP against human serum. Identity has been shown by double diffusion against human serum and a known antibody to human IgG.
Method of Purification	Affinity chromatography
Appearance	Clear colourless solution
Concentration	10,000 U/ml
Volume	1.0 ml
Total Amount	0.75 mg
Buffer	Glycine Buffered Saline pH 7.4

Sterility	0.2µm sterile filtered
Date of manufacture	29-Apr-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 -20 °C

Signed:

Date: 02-Sep-2014



Production Biologist