

Certificate of Analysis

CloneDetect

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

K8215

Batch #	CDD1019A
----------------	-----------------

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 Ruminants; 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
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: 01-NOV-2019

Kieu-Khiem Nguyen

Production Biologist