

Title: **Quality Control Record Sheet**
AQP4 Ab ELISA Version 2 Kit QC Data

Document: wi/qcrs/0.36a
Revision: 2
Date: 8.10.18
Page 1 of 1

Prepared by: G. Flood
Approved by: S. Stephenson

Kit lot number: 2KAQE66

Kit expiry date: 23rd February 2022

Reagent Lot Numbers:

AQP4 Coated Wells: 2AQE72

Expiry date: 11 October 2022

AQP4-Biotin: 2AOB26

Expiry date: 23 November 2022

Reconstitution

Buffer for AQP4-Biotin: 2AOBR22

Expiry date: 03 July 2022

SA-POD: GEP414

Expiry date: 24 August 2022

Diluent for SA-POD: PB164

Expiry date: 01 March 2023

Peroxidase Substrate: PTS126

Expiry date: 23 November 2023

Stop Solution: HESP110

Expiry date: 01 March 2023

Concentrated Wash Solution: AQCW55

Expiry date: 20 October 2022

Calibrators: 2AQEC71

Expiry date: 25 February 2022

Negative Control: 2AQENC71

Expiry date: 25 February 2022

Positive Controls: 2AQEPCa/b71

Expiry date: 25 February 2022

Calibrator Curve and Controls:

Incubation temp: 21°C

Calibrator Concentration (u/mL)	Absorbance at 450 nm
Negative Control	0.047
1.5	0.084
5	0.311
20	1.422
40	2.396
80*	5.447

*Absorbance derived from reading at 405 nm

Positive Control	Absorbance at 450 nm	Concentration (u/mL)	Range (u/mL)
Control I	1.016	12.4	14 ± 5
Control II	1.607	25.2	25 ± 9

Materials of human origin used in the manufacture of this product have been tested and found non-reactive for HIV1 and 2 and HCV antibodies and HBsAg at the time of testing.

Assay performed by: Becky Coles
Signature: *Becky Coles*
Position: Senior Scientist
Date: 04 May 2021

Authorised by: *Dr G Flood*
Signature: *G Flood*
Position: HEAD OF QC
Date: 24th MAY 2021