

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

Document: **wi/qcrs/0.28a**  
Revision: **3**  
Date: **21 September 2018**  
Page **1 of 1**

Prepared by: **G FLOOD**  
Approved by: **A SALVAGE**

Kit lot number: 2KIE179A Kit expiry date: 30 March 2026  
Kit manufacture date: 31 March 2025

**Reagent Lot Numbers:**

IA-2 Coated Wells: <u>IE240</u>	Expiry date: <u>08 JUN 2026</u>
IA-2-Biotin: <u>IEB168</u>	Expiry date: <u>26 MAY 2026</u>
Recon. Buffer for IA-2-Biotin: <u>IBB187</u>	Expiry date: <u>17 JUN 2026</u>
Streptavidin Peroxidase: <u>GEP691</u>	Expiry date: <u>10 AUG 2026</u>
Diluent for SA-POD: <u>PB184</u>	Expiry date: <u>26 NOV 2026</u>
Reaction Enhancer: <u>IES130</u>	Expiry date: <u>03 OCT 2026</u>
Peroxidase Substrate: <u>PTS191</u>	Expiry date: <u>11 NOV 2028</u>
Stop Solution: <u>HESP137</u>	Expiry date: <u>27 NOV 2026</u>
Concentrated Wash Solution: <u>CW423</u>	Expiry date: <u>10 JUN 2027</u>
Calibrators: <u>IC179a</u>	Expiry date: <u>30 MAR 2026</u>
Controls: <u>INC179a / IPC179a</u>	Expiry date: <u>30 MAR 2026</u>

**Curve and Controls:** Incubation temp: +4 °C OVERNIGHT +21 °C

Calibrator Concentration (units per mL)	Absorbance at 450 nm	Absorbance at 405 nm
Negative control	0.028	0.009
7.5	0.119	0.035
35	0.353	0.103
120	1.884	0.545
350	>4	1.841
4000	>4	2.964

Control	Absorbance at 450 nm	Absorbance at 405 nm	Concentration at 450 nm (units per mL)	Concentration at 405 nm (units per mL)	Range (units per mL)
Positive I	1.930	0.557	121.8	121.5	125 ± 40

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: pp Rebecca Coles  
Signature: *Rebecca Coles*  
Position: Principal Scientist  
Date: 18/19 February 2025

Authorised by: DR G FLOOD  
Signature: *G Flood*  
Position: Head of Quality Control  
Date: 31<sup>ST</sup> MAR 2025