

Title: **Quality Control Record Sheet  
GAD Ab ELISA Kit QC Data**

Document: **wi/qcrs/0.27**  
Revision: **11**  
Date: **10/10/14**  
Page **1 of 1**

Prepared by: **G Flood**  
Approved by: **L Carr**

Kit lot number: KGE529 Kit expiry date: 05 January 2026

Kit manufacture date: 06 January 2025

GAD <sub>65</sub> Coated Wells:	Lot No. <u>GE562</u>	Expiry Date <u>20 April 2026</u>
GAD <sub>65</sub> –Biotin :	Lot No. <u>GEB497</u>	Expiry Date <u>17 April 2026</u>
Recon. Buffer for GAD <sub>65</sub> -Biotin:	Lot No. <u>GBB194</u>	Expiry Date <u>13 April 2026</u>
Streptavidin Peroxidase:	Lot No. <u>GEP670</u>	Expiry Date <u>14 April 2026</u>
Diluent for SA-POD:	Lot No. <u>PB183</u>	Expiry Date <u>16 September 2026</u>
Peroxidase Substrate:	Lot No. <u>PTS189G</u>	Expiry Date <u>09 May 2028</u>
Stop Solution:	Lot No. <u>HESP135</u>	Expiry Date <u>23 September 2026</u>
Concentrated Wash Solution:	Lot No. <u>CW413</u>	Expiry Date <u>07 April 2027</u>
Calibrators:	Lot No. <u>GEC526</u>	Expiry Date <u>05 January 2026</u>
Controls:	Lot No. <u>GENC526/GEPC526</u>	Expiry Date <u>05 January 2026</u>

**Calibrator Curve and Controls:**

Incubation temp: +20 °C

Calibrator Concentration (units per mL)	Absorbance at 450 nm	Absorbance at 405 nm
Negative control	0.020	0.008
5	0.174	0.052
18	0.513	0.152
35	0.940	0.280
120	2.470	0.721
250	3.986	1.209
2000	>4	1.940

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	1.216	0.357	47.7	48.3	50 ± 15

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:

Xing

Signature:

Zhang Xing Feng

Position:

Laboratory Technician

Date:

14 November 2024

Authorised by:

DR G Flood

Signature:

[Signature]

Position:

HEAD OF QC

Date:

18<sup>TH</sup> DEC 2024