

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: KGE510A Kit expiry date: 26 May 2025

Kit manufacture date: 27 May 2024

GAD ₆₅ Coated Wells:	Lot No. <u>GE543</u>	Expiry Date <u>04 September 2025</u>
GAD ₆₅ -Biotin :	Lot No. <u>GEB477</u>	Expiry Date <u>03 September 2025</u>
Recon. Buffer for GAD ₆₅ -Biotin:	Lot No. <u>GBB188</u>	Expiry Date <u>06 September 2025</u>
Streptavidin Peroxidase:	Lot No. <u>GEP623</u>	Expiry Date <u>10 September 2025</u>
Diluent for SA-POD:	Lot No. <u>PB180</u>	Expiry Date <u>18 January 2026</u>
Peroxidase Substrate:	Lot No. <u>PTS187E</u>	Expiry Date <u>30 November 2027</u>
Stop Solution:	Lot No. <u>HESP130</u>	Expiry Date <u>08 February 2026</u>
Concentrated Wash Solution:	Lot No. <u>CW386</u>	Expiry Date <u>25 July 2026</u>
Calibrators:	Lot No. <u>GEC507a</u>	Expiry Date <u>26 May 2025</u>
Controls:	Lot No. <u>GENC507a/GEPC507a</u>	Expiry Date <u>26 May 2025</u>

Calibrator Curve and Controls:

Incubation temp: +21 °C

Calibrator Concentration (units per mL)	Absorbance at 450 nm	Absorbance at 405 nm
Negative control	0.014	0.003
5	0.172	0.051
18	0.555	0.163
35	0.911	0.269
120	2.761	0.803
250	>4	1.273
2000	>4	2.021

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	1.315	0.386	50.4	51.5	49 ± 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:

Signature:

Position:

Date:

Xing

Zhang Ringfeng

Laboratory Technician

02 April 2024

Authorised by:

Signature:

Position:

Date:

DR G FLOOD

G Flood

HEAD OF QC

17TH MAY 2024