

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

Kit lot number: **KGE431** Kit expiry date: **22 August 2022**

Kit manufacture date: **23 August 2021**

GAD ₆₅ Coated Wells:	Lot No. <u> GE467 </u>	Expiry Date <u> 27 December 2022 </u>
GAD ₆₅ -Biotin :	Lot No. <u> GEB401 </u>	Expiry Date <u> 24 December 2022 </u>
Recon. Buffer for GAD ₆₅ -Biotin:	Lot No. <u> GBB163 </u>	Expiry Date <u> 21 December 2022 </u>
Streptavidin Peroxidase:	Lot No. <u> GEP434 </u>	Expiry Date <u> 05 January 2023 </u>
Diluent for SA-POD:	Lot No. <u> PB166 </u>	Expiry Date <u> 01 June 2023 </u>
Peroxidase Substrate:	Lot No. <u> PTS175L </u>	Expiry Date <u> 13 April 2023 </u>
Stop Solution:	Lot No. <u> HESP112 </u>	Expiry Date <u> 22 June 2023 </u>
Concentrated Wash Solution:	Lot No. <u> CW289 </u>	Expiry Date <u> 06 January 2024 </u>
Calibrators:	Lot No. <u> GEC432 </u>	Expiry Date <u> 22 August 2022 </u>
Controls:	Lot No. <u> GENC432/GEPC432 </u>	Expiry Date <u> 22 August 2022 </u>

Calibrator Curve and Controls:

Incubation temp: **+23 °C**

Calibrator Concentration (units per mL)	Absorbance at 450 nm	Absorbance at 405 nm
Negative control	0.010	0.003
5	0.196	0.061
18	0.571	0.170
35	0.988	0.299
120	2.611	0.780
250	3.547	1.083
2000	>4	2.042

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	1.346	0.399	48.6	47.9	51 ± 10

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: *Zhong Qiang Feng*
Position: **Laboratory Technician**
Date: **02 August 2021**

Authorised by: **DR G FLOOD**
Signature: *G Flood*
Position: **HEAD OF QC**
Date: **6TH AUG 2021**