

Title: **Quality Control Record Sheet**
RBA AChRab Assay QC data

Document: **wi/qcrs/0.14a**
Revision: **7**
Date: **6 Sep 2018**
Page **1 of 2**

Prepared by: Geoff Flood
Approved by: Rachael Price

Kit lot no:	<u>KRBA317/A</u>	Expiry date:	<u>24 Jan 2022</u>
Labelled Receptor lot no:	<u>RBA317</u>	Expiry date:	<u>24 Jan 2022</u>
Reconstitution buffer lot no:	<u>164B</u>	Expiry date:	<u>16 Apr 2022</u>
Precipitation enhancer lot no:	<u>PE180</u>	Expiry date:	<u>04 Aug 2022</u>
Anti human IgG lot no:	<u>179Gt</u>	Expiry date:	<u>28 Aug 2022</u>
Normal Human serum lot no:	<u>DLA116k</u>	Expiry date:	<u>28 Aug 2022</u>
Washing solution lot no:	<u>411W</u>	Expiry date:	<u>01 Sep 2022</u>
Negative control lot no:	<u>124Naa</u>	Expiry date:	<u>28 Aug 2022</u>
Positive control lot no:	<u>116PCb</u>	Expiry date:	<u>28 Aug 2022</u>

Total cpm in 50 μ L of labelled receptor: 72345

Serum sample (5 μ L)	cpm bound	nmole/litre toxin bound
Negative control	1226	-
Positive control (range)	9688	5.0 (2.8 – 6.4)
QC sera (neat and diluted in normal human serum)		
(A) K3 (lot <u> B </u>)	5350	2.5
K3/2	3943	1.6
(B) K4 (lot <u> E </u>)	2302	0.64
K4/2	1849	0.37
(C) K5 (lot <u> C </u>)	12776	6.9
K5/2	8457	4.3
ϵ -specific serum	7513	3.7
MG7a	17994	10.0
MG6b	14670	8.0

Specific activity of toxin (K): 209 Ci/mmol % Counter Efficiency: 72.3%

Receptor labelling date: 22 Nov 2021 Receptor expiry date: 24 Jan 2022