

Title: **Quality Control Record Sheet**
RBA AChRab Assay QC data
 Prepared by: Geoff Flood
 Approved by: Rachael Price

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

Kit lot no:	<u>KRBA312/A</u>	Expiry date:	<u>06 Sep 2021</u>
Labelled Receptor lot no:	<u>RBA312</u>	Expiry date:	<u>06 Sep 2021</u>
Reconstitution buffer lot no:	<u>163B</u>	Expiry date:	<u>30 Jan 2022</u>
Precipitation enhancer lot no:	<u>PE178</u>	Expiry date:	<u>14 Apr 2022</u>
Anti human IgG lot no:	<u>179Gk</u>	Expiry date:	<u>11 Apr 2022</u>
Normal Human serum lot no:	<u>DLA116d</u>	Expiry date:	<u>11 Apr 2022</u>
Washing solution lot no:	<u>405W</u>	Expiry date:	<u>14 Apr 2022</u>
Negative control lot no:	<u>124Nu</u>	Expiry date:	<u>11 Apr 2022</u>
Positive control lot no:	<u>114PCak</u>	Expiry date:	<u>11 Apr 2022</u>

Total cpm in 50µL of labelled receptor: 77250

Serum sample (5 µL)	cpm bound	nmole/litre toxin bound
Negative control	1207	-
Positive control (range)	9579	5.8 (3.3 – 6.7)
QC sera (neat and diluted in normal human serum)		
(A) K3 (lot <u>B</u>)	4851	2.5
K3/2	3559	1.6
(B) K4 (lot <u>E</u>)	2034	0.57
K4/2	1757	0.38
(C) K5 (lot <u>C</u>)	11544	7.2
K5/2	7894	4.6
ε-specific serum	8354	4.9
MG7a	16126	10.3
MG6b	13340	8.4

Specific activity of toxin (K): 177 Ci/mmol % Counter Efficiency: 73.4%
 Receptor labelling date: 05 Jul 2021 Receptor expiry date: 06 Sep 2021