

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>KRBA307/A</u>	Expiry date:	<u>19 Apr 2021</u>
Labelled Receptor lot no:	<u>RBA307</u>	Expiry date:	<u>19 Apr 2021</u>
Reconstitution buffer lot no:	<u>161B</u>	Expiry date:	<u>26 Aug 2021</u>
Precipitation enhancer lot no:	<u>PE176</u>	Expiry date:	<u>22 Oct 2021</u>
Anti human IgG lot no:	<u>179Gd</u>	Expiry date:	<u>21 Nov 2021</u>
Normal Human serum lot no:	<u>DLA111w</u>	Expiry date:	<u>21 Nov 2021</u>
Washing solution lot no:	<u>400W</u>	Expiry date:	<u>25 Nov 2021</u>
Negative control lot no:	<u>124Nm</u>	Expiry date:	<u>21 Nov 2021</u>
Positive control lot no:	<u>114PCad</u>	Expiry date:	<u>21 Nov 2021</u>

Total cpm in 50 $\mu$ L of labelled receptor: 76076

Serum sample (5 $\mu$ L)	cpm bound	nmole/litre toxin bound
Negative control	1114	-
Positive control (range)	8761	4.9 (3.3 – 6.7)
QC sera (neat and diluted in normal human serum)		
(A) K3 (lot <u>      B      </u> )	4813	2.4
K3/2	3344	1.4
(B) K4 (lot <u>      E      </u> )	2082	0.62
K4/2	1666	0.35
(C) K5 (lot <u>      C      </u> )	11783	6.8
K5/2	7909	4.4
$\epsilon$ -specific serum	7569	4.1
MG7a	16241	9.7
MG6b	14018	8.3

Specific activity of toxin (K): 192 Ci/mmol % Counter Efficiency: 73.4%

Receptor labelling date: 15 Feb 2021 Receptor expiry date: 19 Apr 2021