

Title: **Quality Control Record Sheet
TRAb Coated Tubes QC Data**

Document: **wi/qcrs/0.23**
 Revision: **4**
 Date: **11.08.10**
 Page **1 of 1**

Prepared by: G Flood
 Approved by: W M Williams

Kit Lot Number: KTCT357

Kit expiry date: 10 Mar 2021

125I -TSH Tracer:	Lot No.	TCT321	Expiry Date	10 Mar 2021
TRAb Assay Coated Tubes:	Lot No.	2RC946	Expiry Date	15 May 2021
Calibrators:	Lot No.	2CTC232	Expiry Date	20 Oct 2022
Negative Control:	Lot No.	2CTC232N	Expiry Date	20 Oct 2022
Positive Control:	Lot No.	2CTC232P	Expiry Date	20 Oct 2022
Wash Buffer:	Lot No.	TEW134	Expiry Date	29 Sep 2023
Start Buffer:	Lot No.	2aAD106	Expiry Date	29 Mar 2022

Total counts: 85140 cpm

Temp. 21 °C

	<u>u/Litre</u>	<u>% bound</u>	<u>% inhibition</u>	
Calibrator				
1	1	12.12	14	
2	2	10.40	26	
3	8	5.50	61	
4	40	1.64	88	
CONTROL	% bound	% inhibition	u/Litre	Range
I (Negative)	14.09	-	-	
II (Positive)	8.24	41	3.7	4.2 +/- 1.2 u/L

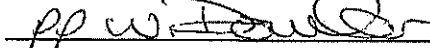
Materials of human origin used in the manufacture of this product have been tested and found non-reactive for HIV1 and 2 and HCV antibodies and HBsAg at the time of testing.

Assay Date: 06 Jan 2021

Authorised by: SARAH NUTE

Performed by: Claire Fisher

Signature: 

Signature: 

Position: pp Head of Quality Control

Position: Laboratory Technician

Date: 12 JAN 2021