

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

Document: w/qcrs/0.23  
 Revision: 5  
 Date: 11.05.21  
 Page: 1 of 1

Prepared by: G Flood  
 Approved by: W Dowler

Kit Lot Number: KTCT375

Kit expiry date: 27 Jul 2022

125I -TSH Tracer:	Lot No.	TCT339	Expiry Date	27 Jul 2022
TSH Receptor Coated Tubes:	Lot No.	2RC990	Expiry Date	14 Sep 2022
Calibrators:	Lot No.	2CTC250	Expiry Date	02 Mar 2024
Negative Control:	Lot No.	2CTC250N	Expiry Date	02 Mar 2024
Positive Control:	Lot No.	2CTC250P	Expiry Date	02 Mar 2024
Concentrated Wash Solution:	Lot No.	TEW138	Expiry Date	06 Dec 2024
Start Buffer:	Lot No.	2aAD108	Expiry Date	19 Jun 2023

Total counts: 85317 cpm

Temp. 21 °C

	<u>IU/L</u>	<u>% bound</u>	<u>% inhibition</u>	
<b>Calibrator</b>				
1	1	13.63	18	
2	2	12.26	26	
3	8	6.24	62	
4	40	1.49	91	
<b>CONTROL</b>	<u>% bound</u>	<u>% inhibition</u>	<u>IU/L</u>	<u>Range</u>
I (Negative)	16.58	-	-	
II (Positive)	9.26	44	4.2	4.2 +/- 1.2 IU/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: 26 May 2022

Authorised by: SARAH NUTE

Performed by: Morwen Williams

Signature: 

Signature: 

Position: Head of Quality Control

Position: Scientist

Date: 27 MAY 2022.