

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

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

Prepared by: G Flood  
Approved by: W Dowler

Kit Lot Number: KTCT363A

Kit expiry date: 25 Aug 2021

125 <sub>I</sub> -TSH Tracer:	Lot No.	TCT327	Expiry Date	25 Aug 2021
TSH Receptor Coated Tubes:	Lot No.	2RC961	Expiry Date	09 Nov 2021
Calibrators:	Lot No.	2CTC238	Expiry Date	23 Mar 2023
Negative Control:	Lot No.	2CTC238N	Expiry Date	23 Mar 2023
Positive Control:	Lot No.	2CTC238P	Expiry Date	23 Mar 2023
Concentrated Wash Solution:	Lot No.	TEW134	Expiry Date	29 Sep 2023
Start Buffer:	Lot No.	2aAD106	Expiry Date	29 Mar 2022

Total counts: 85501 cpm

Temp. 21 °C

	IU/L	% bound	% inhibition	
<b>Calibrator</b>				
1	1	14.46	17	
2	2	12.73	27	
3	8	6.28	64	
4	40	1.68	90	
<b>CONTROL</b>	<b>% bound</b>	<b>% inhibition</b>	<b>IU/L</b>	<b>Range</b>
I (Negative)	17.41	-	-	
II (Positive)	9.47	46	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: 23 Jun 2021

Authorised by: SARAH NUTE

Performed by: Wendy Dowler

Signature: SRNWK

Signature: pp Menni

Position: Head of Quality Control

Position: Section Leader

Date: 30 JUN 2021