

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: KTCT370A

Kit expiry date: 10 Mar 2022

125I -TSH Tracer:	Lot No.	TCT334	Expiry Date	10 Mar 2022
TSH Receptor Coated Tubes:	Lot No.	2RC979	Expiry Date	28 May 2022
Calibrators:	Lot No.	2CTC245	Expiry Date	17 Aug 2023
Negative Control:	Lot No.	2CTC245N	Expiry Date	17 Aug 2023
Positive Control:	Lot No.	2CTC245P	Expiry Date	17 Aug 2023
Concentrated Wash Solution:	Lot No.	TEW136	Expiry Date	18 Apr 2024
Start Buffer:	Lot No.	2aAD107	Expiry Date	18 Oct 2022

Total counts: 85253 cpm

Temp. 21 °C

	IU/L	% bound	% inhibition	
Calibrator				
1	1	11.67	18	
2	2	10.28	28	
3	8	4.98	65	
4	40	1.39	90	
CONTROL	% bound	% inhibition	IU/L	Range
I (Negative)	14.23	-	-	
II (Positive)	7.98	44	3.8	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: 06 Jan 2022

Authorised by: SARAH NUTE

Performed by: Hannah Sanyaolu

Signature: [Signature]

Signature: [Signature]

Position: Head of Quality Control

Position: Senior Scientist

Date: 11 JAN 2022