

Invoice Date & Tax Point 03 February 2021		Invoice No. 33121-21		RSR LIMITED TEL:+44(29)2068 9299 PARC TY GLAS FAX: +44(29)2075 7770 LLANISHEN CARDIFF CF14 5DU VAT REGISTRATION NO: GB 736 8339 02 EORI NO: GB 736 8339 02 000		
Date Requested 14, 22 & 26 January 2021		Customer Purchase No. A12334, 14383, A14384, A15912				
Bill to:- Cisbio Bioassays Parc Marcel Boiteux BP 84175 30 200 Codolet FRANCE			Deliver to:- SDV CDG Roissy Airport For: Cisbio Bioassays BP 32 91192 Gif Sur Yvette Cedex FRANCE			
VAT No.		FR 463 122 618 94		CCN 3822 0000		
Quantity	Order Code	Description	Unit Price EUROS	Total Amount		
1	AQP4/96/2	AQP4 Ab ELISA v2 96W	€ 450.00	450.00		
1	MSK/TRS	MuSK Tracer (inc. Control Set)	€ 145.00	145.00		
1	RBAN/25	AChR Ab IPA 25T	€ 42.00	42.00		
1	LEM/25	VGCC Ab (LEMS) RIA 25T	€ 220.00	220.00		
4	VGK/25	VGKC Ab RIA 25T	€ 227.00	908.00		
Total MBq of 125-I = 0.174						
"I the undersigned do hereby certify that the goods mentioned on the invoice are wholly of United Kingdom origin and that the particulars covered are true and correct." For and on behalf of RSR Limited. (Authorised Signatory) <i>R. Man</i>			On arrival keep refrigerated 2-8°C		Subtotal	1,765.00
					Shipping and Handling	0.00
					Invoice Total	EUROS
Shipping Date		03 February 2021				
Our bankers:		HSBC Bank Plc GB44HBUK40127639544770 HBUKGB4B		Account Name: RSR Limited Sort Code: 40-12-76 Account No.: 39544770		
Shipper's Checker (Intl) GM	Shipping Details		Weight: Net:	3 Kg		
<i>GM</i> Shipper's Signature RH <i>R. Man</i>	Courier: Alexian Flight: LX 0319 04/02/21 LX 0632 06/02/21		Gross:	4 Kg		
			Dimensions (cm):	(Qty)	SBN	
			41 x 30 x 40	(x 1)	3	
		1 pieces				

Fact Re 08/02/2021

M. BONDIS

Title: **Quality Control Record Sheet**
AQP4 Ab ELISA Version 2 Kit QC Data

Document: **wi/qcrs/0.36a**
 Revision: **2**
 Date: **8.10.18**
 Page **1 of 1**

Prepared by: **G. Flood**
 Approved by: **S. Stephenson**

Kit lot number: 2KAQE64

Kit expiry date: 31st October 2021

Reagent Lot Numbers:

AQP4 Coated Wells: 2AQE70

Expiry date: 01 May 2022

AQP4-Biotin: 2AQB24b

Expiry date: 28 April 2022

Reconstitution

Buffer for AQP4-Biotin: 2AQBR22

Expiry date: 03 July 2022

SA-POD: GEP380

Expiry date: 14 January 2022

Diluent for SA-POD: PB161

Expiry date: 18 October 2022

Peroxidase Substrate: PTS124

Expiry date: 03 September 2023

Stop Solution: HESP107

Expiry date: 15 September 2022

Concentrated Wash Solution: AQCW54

Expiry date: 10 July 2022

Calibrators: 2AQEC68

Expiry date: 10 November 2021

Negative Control: 2AQENC68

Expiry date: 10 November 2021

Positive Controls: 2AQEPCa/b68

Expiry date: 10 November 2021

Calibrator Curve and Controls:

Incubation temp: 21°C

Calibrator Concentration (u/mL)	Absorbance at 450 nm
Negative Control	0.042
1.5	0.066
5	0.236
20	1.013
40	2.466
80*	5.307

*Absorbance derived from reading at 405 nm

Positive Control	Absorbance at 450 nm	Concentration (u/mL)	Range (u/mL)
Control I	0.527	11.9	11 ± 4
Control II	1.361	25.2	26 ± 9

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 performed by: Sinead Stephenson

Authorised by: DR G FLOOD

Signature: S. Stephenson

Signature: [Handwritten Signature]

Position: Lab Technician

Position: HEAD OF QC

Date: 15 January 2021

Date: 22ND JAN 2021

Title: **Quality Control Record Sheet
VGCC Ab QC Data**

Prepared by: G Flood
Approved by: J Davies

Document: **wi/qcrs/0.15**
Revision: **8**
Date: **15/08/18**
Page **1 of 1**

Kit Lot Number: LK225 Kit Expiry Date: 09 APRIL 2021

Kit Manufacture Date: 22 JANUARY 2021

¹²⁵I-VGCC for total binding lot number: LT225 Expiry Date: 09 APRIL 2021

¹²⁵I-VGCC for non-specific binding lot number: NS225 Expiry Date: 09 APRIL 2021

Dilution buffer lot number: 219LB Expiry Date: 09 APRIL 2021

Anti-human IgG lot number: 25LGae Expiry Date: 09 APRIL 2021

Wash solution lot number: 142LW Expiry Date: 09 APRIL 2021

Negative control lot number: 19LNh Expiry Date: 09 APRIL 2021

Positive control lot number: 24LPn Expiry Date: 09 APRIL 2021

At Manufacture Date: 6436 dpm = 1 fmole


At QC Assay Date¹: 6146 dpm = 1 fmole (conversion factor A)


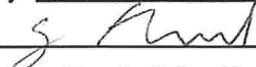
Counter Efficiency: 76.25 % (B) ² Positive control range: 160 ± 70 pmol/L

Sample	NSB		T			
	Mean cpm - background	fmol/ Pellet (E)	Mean cpm - background	fmol/ pellet (D)	¹²⁵ I-conotoxin labelled VGCC bound (pmol/L serum)	pmol/L serum VGCC Ab (after subtraction of negative control)
Total counts	17606	3.757	17990	3.839		
Negative control	243	0.052	459	0.098	18.5(G)	-
Positive control	278	0.059	2260	0.482	169.2 (F)	150.7 ² (H)
Test sample 1	307	0.066	2822	0.602	214.7 (F)	196.2 (H)
Test sample 2b	234	0.050	1111	0.237	74.9 (F)	56.4 (H)
Test sample 3b	286	0.061	2387	0.509	179.4 (F)	160.9 (H)
Test sample 4a	162	0.035	1843	0.393	143.5 (F)	125.0 (H)
Test sample 5d	261	0.056	1823	0.389	133.4 (F)	114.9 (H)

Assay date (weeks after QC assay)	Decay factor (C)	Assay date (weeks after QC assay)	Decay factor (C)
Up to 1 week	1.00	+ 6 – 7 weeks	0.61
+ 1 – 2 weeks	0.92	+ 7 – 8 weeks	0.56
+ 2 – 3 weeks	0.85	+ 8 – 9 weeks	0.52
+ 3 – 4 weeks	0.78	+ 9 – 10 weeks	0.48
+ 4 – 5 weeks	0.72	+ 10 – 11 weeks	0.44
+ 5 – 6 weeks	0.66		

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.

QC Assay Date¹: 26 JANUARY 2021
Performed by: L GUO
Signature: 
Position: Senior Scientist

Authorised by: 
Signature: 
Position: Head of Quality Control
Date: 29th JAN 2021

Title: **Quality Control Record Sheet
VGKC Ab Kit QC Data**

Document: **wi/qcrs/0.37**
Revision: **3**
Date: **15.08.18**
Page **1 of 1**

Prepared by: G Flood
Approved by: S Stephenson

Kit lot number: VGK76 Kit Expiry Date: 02 Apr 2021
¹²⁵I-Labelled VGKC lot number: VKT76 Expiry Date: 02 Apr 2021
 Anti-Human IgG lot number: VKG12 Expiry Date: 15 Jan 2022
 Assay Buffer lot number: VKB53 Expiry Date: 05 Jul 2021
 Negative Control: VKNC73 Expiry Date: 30 Apr 2021
 Positive Controls I and II: VKPCa/b73 Expiry Date: 30 Apr 2021

Tracer manufacture date: 22nd January 2021
 Total cpm in 50 µL of ¹²⁵I – VGKC: 54792
 Specific activity of toxin (K): 3463 Ci/mmol
 Counter efficiency: 72.3 %

Assay date (weeks after manufacture)	Decay factor (A)
Up to 1 week	1.0
+1 – 2 weeks	1.1
+2 – 3 weeks	1.2
+3 - 4 weeks	1.3
+4 – 5 weeks	1.4
+5 – 6 weeks	1.5
+6 – 7 weeks	1.6
+7 – 8 weeks	1.75
+8 – 9 weeks	1.9
+9 – 10 weeks	2.1

	Mean cpm	pM	Range (pM)
Negative Control	1043	-	-
Positive Control I	12641	442.0	450 ± 130
Positive Control II	6661	214.1	210 ± 60

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 performed by: Fiona Ryrie
 Signature: Fiona Ryrie
 Position: Scientific Assistant
 Date of assay: 27th January 2021

Authorised by: DR G Flood
 Signature: G Flood
 Position: HEAD OF QC
 Date: 27th JAN 2021