



# PRODUCT LIST 2019



Immunoassays  
RIA & ELISA



## AUTO-IMMUNITY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (months)
<b>ACHRAB ELISA</b>	Microplate ACE/96	ELISA	96	Serum 100 µL	0.5-20 nmol/mL	16-20h 2-8°C + 1h + 1h + 30 min RT Shaking	9
<b>AQP4 AB ELISA</b>	Microplate AQP4/96	ELISA	96	Serum 50 µL	5-160 U/mL	2h + 20 min RT Shaking	9
<b>GADAB ELISA</b>	Microplate GADAB-ELISA	ELISA	96	Serum 25 µL	5-2000 U/mL	1+1h RT Shaking	12
<b>IA-2 AB ELISA</b>	Microplate IA2AB-ELISA	ELISA	96	Serum 50µL	7.5-4000 U/mL	16-20h + 1h 2-8°C +20 min RT Shaking	12
<b>21-OH AB ELISA</b>	Microplate 21E/96	ELISA	96	Serum 50 µL	0.3-100 U/mL	16-20h 2-8°C + 1h + 20 min RT Shaking	6
<b>2 SCREEN ELISA</b>	Microplate 2GI/96	ELISA	96	Serum 50 µL	4-450 U/mL	16-20h 2-8°C + 1h + 20 min RT Shaking	12
<b>TGAB ELISA</b>	Microplate TAE/96/2	ELISA	96	Serum 50 µL	65-5000 U/mL	15 + 15 min RT Shaking	12
<b>TPOAB ELISA</b>	Microplate PE/96/2	ELISA	96	Serum 50 µL	10-5000 U/mL	15 + 15 min RT Shaking	12
<b>ZNT8 AB ELISA</b>	Microplate ZnT8/96	ELISA	96	Serum 25 µL	10-2000 U/mL	16-20h + 1h 2-8°C + 20 min RT Shaking	6

## ONCOLOGY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (months)
<b>CHROMOA</b>	Microplate CGA-ELISA	ELISA	96	Serum/Plasma 20 µL	7-820 ng/mL	16-20h RT + 2h RT Shaking ou 2+2h RT Shaking	20

## OTHER PRODUCTS

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (months)
<b>HYALURONAN</b>	Microplate HYAL-ELISA	ELISA	96	Serum/Plasma 10 µL	2.7-1000 ng/mL	2h + 30 min RT Shaking	12
<b>HYALURONAN</b>	Microplate HYAL-US *	ELISA	96	Serum/Plasma 10 µL	2.7-1000 ng/mL	2h + 30 min RT Shaking	12
<b>FECAL INFLAMARK (S100 A12)</b>	Microplate F-INFL-EL-US*	ELISA	96	Stool 1 - 5 g	0.21-12 µg/g	30 min + 30 min RT Shaking	14
<b>SERUM INFLAMARK** (S100 A12)</b>	Microplate S-INFL-ELISA*	ELISA	96	Serum 30 µL	21-650 ng/mL	16-20h RT + 2h RT Shaking	14
<b>N TERMINAL PROCOLLAGEN III PEPTIDE</b>	Microplate P3NP-ELISA	ELISA	96	Serum/Plasma 30 µL	0.036-30 µg/L	3h RT Shaking	12
<b>N TERMINAL PROCOLLAGEN III PEPTIDE</b>	Microplate P3NP-EL-US*	ELISA	96	Serum/Plasma 30 µL <sup>(1)</sup>	0.036-30 µg/L	3h RT Shaking	12

\* For Research Use Only, not for use in diagnostic procedures.

\*\* On request

<sup>(1)</sup> Urine possible

## AUTO-IMMUNITY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>ACETYLCHOLINE RECEPTOR AUTOANTIBODIES</b>	Radioreceptor ACHR-AB	RRA	25	Serum/Plasma 5 µL	-	2h RT + 2h or 18-24h 2-8°C	6
<b>ANTI-TSH RECEPTOR AUTOANTIBODIES</b>	Coated tubes TR-AB-CT	RIA	60	Serum 100 µL	0.33-40 U/L	2 + 1h RT Shaking	6
<b>INSULIN AUTOANTIBODIES</b>	PEG AAI	RIP	100	Serum/Plasma 50 µL	-	2h + 10 min RT	6
<b>MUSK AUTOANTIBODIES</b>	Precipitation MSK/TR	RIP	25	Serum 5 µL	-	18h RT + 2h 2-8°C	6
<b>THYROGLOBULIN AUTOANTIBODIES</b>	Coated tubes TGAB I STEP	RIA	100	Serum 20 µL	6-2000 IU/mL	1h30 RT Shaking	8
<b>TPO AUTOANTIBODIES</b>	Coated tubes TPO-AB-CT	RIA	50	Serum 50 µL	8-10000 U/mL	1h RT Shaking	10
<b>VGCC AUTOANTIBODIES</b>	Precipitation (2nd autoantibody) VGC/25	RIA	25	Serum 25 µL	-	1h+1h RT	8
<b>VGKC AUTOANTIBODIES</b>	Precipitation (2nd autoantibody) VGK/25	RIP	25	Serum 50 µL	-	16-20h +1h30 RT	8

## BONE/GROWTH

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>HGH</b>	Coated tubes HGH-RIACT	IRMA	100	Serum 50 µL	0.01-25.2 ng/mL	2h RT Shaking	7
<b>IGF1</b>	Coated tubes IGF1-RIACT	IRMA	100	Serum/Plasma 50 µL	1-1000 ng/mL	2+2h RT Shaking	10
<b>OSTEOCALCIN</b>	Coated tubes OSTEO-RIACT	IRMA	100	Serum/Plasma 50 µL	0.4-300 ng/mL	2h RT Shaking	7
<b>PTH (1)</b>	Solid phase-ELSA system ELSA-PTH	IRMA	96	Serum/Plasma 200 µL	0.7-1500 pg/mL	16-20h RT	6

(1) Quick optional peroperative protocol

## ENDOCRINOLOGY / HYPERTENSION

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>ALDOSTERONE</b>	Coated tubes ALDO-RIACT	RIA	100	Serum/Plasma 200 µL (1)	7-1500 pg/mL	3h RT Shaking	9
<b>ANGIOTENSINE I</b>	Coated tubes REN-CT2	RIA	100	Plasma 500 µL	0.15-25 ng/mL	90 min RT Shaking	7
<b>RENIN</b>	Coated tubes RENINE	IRMA	100	Plasma 300 µL	1-320 pg/mL	3h RT Shaking	6

(1) Urine possible

## DIABETES

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>ANTI-INSULIN AUTOANTIBODIES</b>	PEG AAI	RIP	100	Serum/Plasma 50 µL	-	2h + 10 min RT	6
<b>C-PEPTIDE</b>	Coated tubes IRMA-CPEP	IRMA	100	Serum/Plasma (1) 50 µL	0.012-6 nmol/L	2h RT Shaking	8
<b>INSULIN</b>	Coated tubes BI-INS-IRMA	IRMA	100	Serum/Plasma 50 µL	0.2-500 µIU/mL	2h RT Shaking	6

(1) Urine possible

## FERTILITY / REPRODUCTION / HYPOTHALAMIC-PITUITARY AXIS

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>ACTH</b>	Solid phase-ELSA system ELSA-ACTH	IRMA	96	Plasma 200 µL	2-2000 pg/mL	18-22h RT	7
<b>ANDROSTENEDIONE</b>	Coated tubes R-GM-100	RIA	100	Serum 25µL	0.05-10 ng/mL	1h 37°C	9
<b>CORTISOL</b>	Coated tubes CORT-CT2	RIA	100	Serum/Saliva 20 µL /150 µL	20-2000 nmol/L (serum) 1-100 nmol/L (saliva)	2h 37°C (serum) 30 min (saliva)	9
<b>ESTRADIOL</b>	Coated tubes ESTR-CTRIA	RIA	100	Serum/Plasma 100 µL	4-5000 pg/mL	2h30 RT Shaking	9
<b>ESTRADIOL (ULTRASENSITIVE) (1)</b>	Coated tubes ESTR-US-CT	RIA	100	Serum/Plasma 200 µL	5-2000 pmol/L	1h + 1h 37°C	7
<b>HCG</b>	Coated tubesRIAgnost OCPM03-HCG	IRMA	100	Serum/Plasma 100 µL	0-1000 mIU/mL	<i>Qualitative assay</i> 10min + 10min RT Shaking <i>Quantitative assay</i> 30min + 60min RT Shaking	10
<b>FβHCG (Free β chains)</b>	Solid phase-ELSA system ELSA-FβHCG	IRMA	96	Serum/Plasma 100 µL	0.04-50 ng/mL	1+2h RT Shaking	7
<b>17-OH-PROGESTERONE</b>	Coated tubes OHP-CT	RIA	100	Serum/Plasma 25 µL	0.09-75 nmol/L	2h30 37°C	7
<b>PROGESTERONE</b>	Coated tubes PROGEST-RIA	RIA	96	Serum 50 µL	0.05-36 ng/mL	2h 37°C	8
<b>SHBG</b>	Coated tubes SHBG-RIACT	IRMA	100	Serum/Plasma 10 µL	0.5-200 nmol/L	30+30 min RT Shaking	11
<b>TESTOSTERONE</b>	Coated tubes TESTO-CT2	RIA	100	Serum 25 µL	0.3-50 nmol/L	1h 37°C	13

(1) Sensibility 5 pmol/L

## ONCOLOGY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>AFP</b>	Solid phase-ELSA system ELSA2-AFP	IRMA	96	Serum / Plasma / Amniotic fluid 50 µL	0.5-800 ng/mL	30+30 min RT Shaking	7
	Coated tubes RIAGnost OCFA07-AFP	IRMA	100	Serum/Plasma 50 µL	0.3-800 IU/mL	15+15 min RT Shaking	10
<b>CALCITONIN</b>	Coated tubes IRMA-hCT	IRMA	100	Serum 200 µL	1.5-1500 pg/mL	20-24h RT	4
<b>CA 125</b>	Solid phase-ELSA system ELSA-CA 125 II	IRMA	96	Serum/Plasma 100 µL	0.5-500 U/mL	18h RT	7
<b>CA 15-3</b>	Solid phase-ELSA system ELSA-CA 15-3	IRMA	96	Serum/Plasma 20 µL	0.2-240 U/mL	1+1h RT Shaking	7
<b>CA 19-9</b>	Solid phase-ELSA system ELSA-CA 19-9	IRMA	96	Serum/Plasma 100 µL	1.5-240 U/mL	3h 37°C + 3h ou 16-24h RT Shaking	7
<b>CA 50</b>	Coated tubes RIAGnost OCFM07 - CA50	IRMA	100	Serum/Plasma 50 µL	0.4-180 U/mL	2+2h RT Shaking	7
<b>CA 72-4</b>	Solid phase-ELSA system ELSA-CA 72-4	IRMA	96	Serum/Plasma 100 µL	0.8-100 U/mL	4h RT Shaking + 16h-20h 2-6°C	7
<b>CEA</b>	Coated tubes CEA-RIACT	IRMA	100	Serum/plasma 50 µL	1-200 ng/mL	1h RT Shaking	6
<b>CHROMOGRANIN A</b>	Coated tubes CGA-RIACT	IRMA	100	Serum/Plasma 50 µL	1.5-1200 ng/mL	18-20h RT + 2h RT Shaking	9
<b>CYFRA 21-1</b>	Solid phase-ELSA system ELSA-CYFRA-	IRMA	48	Serum 100 µL	0.05-63 ng/mL	18-22h 2-8°C	7
<b>ESTRADIOL (ULTRASENSITIVE) (1)</b>	Coated tubes ESTR-US-CT	RIA	100	Serum/Plasma 200 µL	5-2000 pmol/L	1h + 1h 37°C	7
<b>FBHCG (FREE B CHAINS)</b>	Solid phase-ELSA system ELSA-FBHCG	IRMA	96	Serum/Plasma 100 µL	0.04-50 ng/mL	1+2h RT Shaking	7
<b>HCG</b>	Coated tubesRIAGnost OCPM03-HCG	IRMA	100	Serum/Plasma 100 µL	0-1000 mIU/mL	<i>Qualitative assay</i> 10min + 10min RT Shaking <i>Quantitative assay</i> 30min + 60min RT Shaking	10
<b>NSE</b>	Solid phase-ELSA system ELSA-NSE	IRMA	96	Serum 50 µL	0.3-200 ng/mL	3h RT Shaking	6
<b>PSA</b>	Coated tubes PSA-RIACT	IRMA	100	Serum/Plasma 50 µL	0.3-100 ng/mL	2h RT Shaking	5
<b>PSA (FREE)</b>	Coated tubes FPSA-RIACT	IRMA	50	Serum/Plasma 100 µL	0.16-20 ng/mL	16-20h RT Shaking	5
<b>THYROGLOBULIN</b>	Coated tubes THYRO	IRMA	100	Serum/Plasma 100 µL	0.2-500 ng/mL	3h RT Shaking + 16-20h RT ou 16h RT	6

(1) Sensibility 5 pmol/L

## THYROID

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>THYROGLOBULIN AUTOANTIBODIES</b>	Coated tubes TGAB I STEP	RIA	100	Serum 20 µL	6-2000 IU/mL	1h30 RT Shaking	8
<b>TSH RECEPTOR AUTOANTIBODIES</b>	Coated tubes TR-AB-CT	RIA	60	Serum 100 µL	0.3-40 U/L	2+1h RT Shaking	6
<b>THYRO PEROXYDASE AUTOANTIBODIES</b>	Coated tubes TPO-AB-CT	RIA	50	Serum 50 µL	8-10000 U/mL	1h RT Shaking	10
<b>CALCITONIN</b>	Coated tubes IRMA-hCT	IRMA	100	Serum 200 µL	1.5-1500 pg/mL	20-24h RT	4
<b>T3</b>	Coated tubes RIAGnost OCFE07-T3 / OCFE57-T3	RIA	100 / 500	Serum/Plasma 50 µL	0.1-7.5 ng/mL	45 min RT Shaking	10
Adaptable on RIA-MAT SR300 analyser, for protocol contact us at iva@cisbio.com							
<b>FREE T3</b>	Coated tubes RIAGnost OCFH07-FT3 / OCFH57-FT3	RIA	100 / 500	Serum 50 µL	0.6-50 pg/mL	2h RT Shaking	9
Adaptable on RIA-MAT SR300 analyser, for protocol contact us at iva@cisbio.com							
<b>T4</b>	Coated tubes RIAGnost OCPG07-T4 / OCPG57-T4	RIA	100 / 500	Serum/Plasma 20 µL	2.5-360 ng/mL	45 min RT Shaking	10
Adaptable on RIA-MAT SR300 analyser, for protocol contact us at iva@cisbio.com							
<b>FREE T4</b>	Coated tubes RIAGnost OCFD03-FT4 / OCFD53-FT4	RIA	100 / 500	Serum 100 µL	0.5-70 pg/mL	30+60 min RT Shaking	9
<b>FREE T4</b>	Coated tubes RIAGnost OCFD03-FT4-K / OCFD53-FT4-K	RIA	100 / 500	Serum 50 µL	0.73-70 pg/mL	45 min 37°C Shaking	9
Adaptable on RIA-MAT SR300 analyser, for protocol contact us at iva@cisbio.com							
<b>TBG</b>	PEG RIAGnost OCRU03-TBG	RIA	100	Serum/Plasma 20 µL	0-80 µg/mL	30 min RT	13
<b>THYROGLOBULIN</b>	Coated tubes THYRO	IRMA	100	Serum/Plasma 100 µL	0.2-500 ng/mL	3h RT Shaking + 16-20h RT ou 16-20h RT	6
<b>TSH (1)</b>	Coated tubes RIAGnost OCPL07-TSH / OCPL57-TSH	IRMA	100 / 500	Serum/Plasma 200 µL	0.03-65 µIU/mL	2h RT Shaking	8
Adaptable on RIA-MAT SR300 analyser, for protocol contact us at iva@cisbio.com							

(1) Optional calibrator 100 µIU/mL

## OTHER PRODUCTS

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
<b>GASTRIN</b>	Precipitation (2nd autoantibody) GASK-PR	RIP	100	Serum/Plasma 100 µL	10-1500 µU/mL	2h + 15 min RT	7
<b>N TERMINAL PROCOLLAGEN III PEPTIDE</b>	Coated tubes OCFK07-PIIP	RIA	100	Serum/Plasma 20 µL	0.1-14 U/mL	2+3h RT Shaking	6
<b>TRYPsin</b>	Double antibody immuno- precipitating reagent OCFE07-TRYPS	RIP	100	Serum 100 µL	8-1300 ng/mL	3h + 30 min RT	8

For more information, please visit our website [www.cisbio.com](http://www.cisbio.com)