



PRODUCT LIST 2022



Immunoassays
RIA & ELISA



AUTO-IMMUNITY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (months)
ACHRAB ELISA	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
AQP4 AB ELISA	Microplate AQP4/96	ELISA	96	Serum 50 µL	5-160 U/mL	2h + 20 min RT Shaking	9
GADAB ELISA	Microplate GADAB-ELISA	ELISA	96	Serum 25 µL	5-2000 U/mL	1h + 1h RT Shaking	12
IA-2 AB ELISA	Microplate IA2AB-ELISA	ELISA	96	Serum 50 µL	7.5-4000 U/mL	16-20h + 1h 2-8°C +20 min RT Shaking	12
21-OH AB ELISA	Microplate 21E/96	ELISA	96	Serum 50 µL	0.3-100 U/mL	16-20h 2-8°C + 1h + 20 min RT Shaking	6
2 SCREEN ELISA	Microplate 2GI/96	ELISA	96	Serum 50 µL	4-450 U/mL	16-20h 2-8°C + 1h + 20 min RT Shaking	12
ZNT8 AB ELISA	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)
CHROMOA	Microplate CGA-ELISA	ELISA	96	Serum/Plasma 20 µL	7-820 ng/mL	2h + 2h RT Shaking	20
CHROMOGRANINE A	CGA-ELISA-NG	ELISA	96	Serum/Plasma 20 µL	30.6-1000 ng/mL	1h + 2h RT Shaking	20

FIBROSIS

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (months)
N TERMINAL PROCOLLAGEN III PEPTIDE	Microplate P3NP-ELISA	ELISA	96	Serum/Plasma 30 µL	0.036-30 µg/L	3h RT Shaking	24
N TERMINAL PROCOLLAGEN III PEPTIDE	Microplate P3NP-EL-US*	ELISA	96	Serum/Plasma 30 µL ⁽¹⁾	0.036-30 µg/L	3h RT Shaking	24

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

⁽¹⁾ Urine possible

AUTO-IMMUNITY

	Separation Reference	Method	Number of tests	Sample	Measuring range	Incubation	Shelf life (weeks)
ACETYLCHOLINE RECEPTOR AUTOANTIBODIES	Radioreceptor ACHR-AB	RRA	25	Serum/Plasma 5 µL	-	2h RT + 2h or 18-24h 2-8°C	6
ANTI-TSH RECEPTOR AUTOANTIBODIES	Coated tubes TR-AB-CT	RIA	60	Serum 100 µL	0.33-40 U/L	2 + 1h RT Shaking	6
INSULIN AUTOANTIBODIES	PEG AAI	RIP	100	Serum/Plasma 50 µL	-	2h + 10 min RT	6
MUSK AUTOANTIBODIES	Precipitation MSK/TR	RIP	25	Serum 5 µL	-	18h RT + 2h 2-8°C	6
THYROGLOBULIN AUTOANTIBODIES	Coated tubes TGAB I STEP	RIA	96	Serum 20 µL	6-2000 IU/mL	1h30 RT Shaking	8
TPO AUTOANTIBODIES	Coated tubes TPO-AB-CT	RIA	50	Serum 50 µL	8-10000 U/mL	1h RT Shaking	10
VGCC AUTOANTIBODIES	Precipitation (2nd autoantibody) VGC/25	RIA	25	Serum 25 µL	-	1h+1h RT	8
VGKC AUTOANTIBODIES	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)
HGH	Coated tubes HGH-RIACT	IRMA	100	Serum 50 µL	0.01-25.2 ng/mL	2h RT Shaking	7
IGF1	Coated tubes IGF1-RIACT	IRMA	100	Serum/Plasma 50 µL	1-1000 ng/mL	2+2h RT Shaking	10
OSTEOCALCIN	Coated tubes OSTEO-RIACT	IRMA	100	Serum/Plasma 50 µL	0.4-300 ng/mL	2h RT Shaking	7
PTH (1)	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)
ALDOSTERONE	Coated tubes ALDO-RIACT	RIA	100	Serum/Plasma 200 µL (1)	7-1500 pg/mL	3h RT Shaking	9
ANGIOTENSINE I	Coated tubes REN-CTZP	RIA	96	Plasma 500 µL	0.15-25 ng/mL	90 min RT Shaking	7
RENIN	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)
ANTI-INSULIN AUTOANTIBODIES	PEG AAI	RIP	100	Serum/Plasma 50 µL	-	2h + 10 min RT	6
C-PEPTIDE	Coated tubes IRMA-CPEP	IRMA	100	Serum/Plasma (1) 50 µL	0.012-6 nmol/L	2h RT Shaking	8
INSULIN	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)
ACTH	Solid phase-ELSA system ELSA-ACTH	IRMA	96	Plasma 200 µL	2-2000 pg/mL	18-22h RT	7
ANDROSTENEDIONE	Coated tubes R-GM-100	RIA	100	Serum 25µL	0.05-10 ng/mL	1h 37°C	9
CORTISOL	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
ESTRADIOL	Coated tubes ESTR-CTRIA	RIA	100	Serum/Plasma 100 µL	4-5000 pg/mL	2h30 RT Shaking	9
ESTRADIOL (ULTRASENSITIVE) (1)	Coated tubes ESTR-US-CT	RIA	100	Serum/Plasma 200 µL	5-2000 pmol/L	1h + 1h 37°C	7
HCG	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
FβHCG (Free β chains)	Solid phase-ELSA system ELSA-FβHCG	IRMA	96	Serum/Plasma 100 µL	0.04-50 ng/mL	1+2h RT Shaking	7
17-OH-PROGESTERONE	Coated tubes OHP-CT	RIA	100	Serum/Plasma 25 µL	0.09-75 nmol/L	2h30 37°C	7
PROGESTERONE	Coated tubes PROGEST-RIA	RIA	96	Serum 50 µL	0.05-36 ng/mL	2h 37°C	8
SHBG	Coated tubes SHBG-RIACT	IRMA	100	Serum/Plasma 10 µL	0.5-200 nmol/L	30+30 min RT Shaking	11
TESTOSTERONE	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)
AFP	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
CALCITONIN	Coated tubes IRMA-hCT	IRMA	100	Serum 200 µL	1.5-1500 pg/mL	20-24h RT	4
CA 125	Solid phase-ELSA system ELSA-CA 125 II	IRMA	96	Serum/Plasma 100 µL	0.5-500 U/mL	18h RT	7
CA 15-3	Solid phase-ELSA system ELSA-CA 15-3	IRMA	96	Serum/Plasma 20 µL	0.2-240 U/mL	1+1h RT Shaking	7
CA 19-9	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
CA 50	Coated tubes RIagnost OCFM07 - CA50	IRMA	100	Serum/Plasma 50 µL	0.4-180 U/mL	2+2h RT Shaking	7
CA 72-4	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
CEA	Coated tubes CEA-RIACT	IRMA	100	Serum/plasma 50 µL	1-200 ng/mL	1h RT Shaking	6
CHROMOGRANIN A	Coated tubes CGA-RIACT	IRMA	100	Serum/Plasma 50 µL	1.5-1200 ng/mL	18-20h RT + 2h RT Shaking	9
CYFRA 21-1	Solid phase-ELSA system ELSA-CYFRA-	IRMA	48	Serum 100 µL	0.05-63 ng/mL	18-22h 2-8°C	7
ESTRADIOL (ULTRASENSITIVE) (1)	Coated tubes ESTR-US-CT	RIA	100	Serum/Plasma 200 µL	5-2000 pmol/L	1h + 1h 37°C	7
FBHCG (FREE B CHAINS)	Solid phase-ELSA system ELSA-FBHCG	IRMA	96	Serum/Plasma 100 µL	0.04-50 ng/mL	1+2h RT Shaking	7
HCG	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
NSE	Solid phase-ELSA system ELSA-NSE	IRMA	96	Serum 50 µL	0.3-200 ng/mL	3h RT Shaking	6
PSA	Coated tubes PSA-RIACT	IRMA	100	Serum/Plasma 50 µL	0.3-100 ng/mL	2h RT Shaking	5
PSA (FREE)	Coated tubes FPSA-RIACT	IRMA	50	Serum/Plasma 100 µL	0.16-20 ng/mL	16-20h RT Shaking	5
THYROGLOBULIN	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)
THYROGLOBULIN AUTOANTIBODIES	Coated tubes TGAB I STEP	RIA	96	Serum 20 µL	6-2000 IU/mL	1h30 RT Shaking	8
TSH RECEPTOR AUTOANTIBODIES	Coated tubes TR-AB-CT	RIA	60	Serum 100 µL	0.3-40 U/L	2+1h RT Shaking	6
THYRO PEROXYDASE AUTOANTIBODIES	Coated tubes TPO-AB-CT	RIA	50	Serum 50 µL	8-10000 U/mL	1h RT Shaking	10
CALCITONIN	Coated tubes IRMA-hCT	IRMA	100	Serum 200 µL	1.5-1500 pg/mL	20-24h RT	4
T3	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							
FREE T3	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							
T4	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							
FREE T4	Coated tubes RIAGnost OCFD03-FT4 / OCFD53-FT4	RIA	100 / 500	Serum 100 µL	0.5-70 pg/mL	30+60 min RT Shaking	9
FREE T4	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							
TBG	PEG RIAGnost OCRU03-TBG	RIA	100	Serum/Plasma 20 µL	0-80 µg/mL	30 min RT	13
THYROGLOBULIN	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
TSH (1)	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)
GASTRIN	Precipitation (2nd autoantibody) GASK-PR	RIP	100	Serum/Plasma 100 µL	10-1500 µU/mL	2h + 15 min RT	7
N TERMINAL PROCOLLAGEN III PEPTIDE	Coated tubes OCFK07-PIIP	RIA	100	Serum/Plasma 20 µL	0.1-14 U/mL	2+3h RT Shaking	6
TRYPsin	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