

A close-up photograph of laboratory glassware, including a test tube and a beaker, with a green banner overlaid in the top left corner. The banner contains the text 'NEW 2022 SCHEDULE'.

**NEW**  
2022 SCHEDULE

# Availability Schedule 2022

In-Vitro RIA Products

# 2022 - AVAILABILITY SCHEDULE by product

## In-Vitro RIA Products

### BANK HOLIDAYS IN 2022:

1<sup>st</sup> January, 18<sup>th</sup> April, 1<sup>st</sup> May, 8<sup>th</sup> May, 26<sup>th</sup> May, 6<sup>th</sup> June, 14<sup>th</sup> July,  
15<sup>th</sup> August, 1<sup>st</sup> November, 11<sup>th</sup> November, 25<sup>th</sup> December.  
Please order one day before normal ordering day

**ORDERS:** Phone: +33 169 857 253 or +33 169 857 546  
Fax: +33 169 857 521

PARAMETER	KIT REFERENCE	SHELF-LIFE (weeks)	ACT (kBq)	JANUARY			FEBRUARY			MARCH			APRIL			MAY			JUNE		
				1-9	10-19	20-31	1-9	10-19	20-27	1-9	10-19	20-31	1-9	10-19	20-30	1-9	10-18	19-31	1-9	10-19	20-30
ACTH	ELSA-ACTH	7 weeks	592		We 19			We 16			We 16		We 13		We 11	Mo 30					
AFP	ELSA2-AFP	7 weeks	277,5		We 26			We 23		We 23			Th 21		We 18						
	OCFA07-AFP	10 weeks	300	We 05					We 02				We 27								
ALDOSTERONE	ALDO-RIACT	9 weeks	111		Mo 17					Mo 14				Mo 09							
ANDROSTENEDIONE	R-GM-100	9 weeks	115		We 12		We 09		We 09			We 06		We 04							
ANGIOTENSIN I	REN-CT2P	7 weeks	150			Th 27			Th 24		Th 24			Fr 22		Th 19			Th 16		
ANTI-ACETYLCHOLINE RECEPTOR AUTOANTIBODIES	ACHR-AB	6 weeks	40		We 12		We 02		We 02		We 30		We 27			We 25				We 22	
ANTI-INSULIN ANTIBODIES	AAI	6 weeks	52	Mo 03		Mo 31		Mo 28		Mo 28			Mo 25		Mo 23						
ANTI-THYROGLOBULINE ANTIBODIES	TGAB I STEP	8 weeks	137		Th 13			Th 10		Th 10		Th 07		Th 05			Th 02			Th 30	
ANTI-TPO AUTOANTIBODIES	TPO-AB-CT	10 weeks	40				We 02				We 30				We 25						
ANTI-TSH RECEPTOR AUTOANTIBODIES	TR-AB-CT	6 weeks	100			We 26		We 23		We 23		We 20		We 18			We 15				
ANTI-VGKC AUTOANTIBODIES	VGK/25	8 weeks	30			We 26			We 09					We 04						We 29	
ANTI-VGCC AUTOANTIBODIES	VGC/25	9 weeks	12			We 26			We 09					We 04						We 29	
CALCITONIN	IRMA-HCT	4 weeks	611	We 05	We 19		We 02	We 16	We 02	We 16	We 30		We 13	We 27	We 11	Mo 30					
CA 125	ELSA-CA125II	7 weeks	481	We 05			We 02		We 02		We 30		We 27		We 25						
CA 15-3	ELSA-CA15-3	7 weeks	555			We 26		We 23			We 23		Th 21		We 18						
CA 19-9	ELSA-CA19-9	7 weeks	370			We 26		We 23			We 23		Th 21		We 18						
CA 50	OCFM07-CA50	7 weeks	300	We 05				We 16			We 30				We 11						
CA72-4	ELSA-CA72-4	7 weeks	370			We 26		We 23			We 23		Th 21		We 18						
CEA	CEA-RIACT	6 weeks	350	Mo 03		Mo 31		Mo 28			Mo 28		Mo 25			Mo 23					
CHROMOGRANIN A	CGA-RIACT	9 weeks	407	We 05					We 02				We 27								
CORTISOL	CORT-CT2	9 weeks	250		We 12		We 09		We 09			We 06		We 04		Mo 30					
C-PEPTIDE	IRMA-CPEP	8 weeks	460	We 05				We 16			We 30				We 11						
CYFRA-21-1	ELSA-CYFRA-	7 weeks	277,5			Fr 28			Fr 25		Fr 25			Fr 22		Fr 20					
ESTRADIOL	ESTR-CTR1A	9 weeks	93					Mo 14					Mo 11		We 11						
	ESTR-US-CT	7 weeks	85		Fr 14			Fr 11		Fr 11		Fr 08		Fr 06		Mo 30					
FT3/FT4	OCFH07-FT3	9 weeks	150	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
	OCFH57-FT3	9 weeks	750	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
	OCFD03-FT4	9 weeks	150	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
	OCFD03-FT4-K	9 weeks	150	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
	OCFD53-FT4	9 weeks	750	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
GASTRIN	GASK-PR	7 weeks	37		We 12		We 09		We 09			We 06		We 04		We 25					
HCG - Free β chains	ELSA-FBHCG	7 weeks	555			We 26		We 23		We 23			Th 21		We 18						
HCG	OCPM03-HCG	10 weeks	300				We 16					We 13		We 11							
HGH	HGH-RIACT	7 weeks	308		We 19		We 16			We 16		We 13		We 11							
IGF1	IGF1-RIACT	10 weeks	370		Mo 17						Mo 14			Mo 09							
INSULIN	BI-INS-IRMA	6 weeks	460			Fr 28			Fr 25		Fr 25		Fr 22		Fr 20						
MuSK	MSK/TR	6 weeks	50			We 26		We 23		We 23		We 20		We 18			We 15				
NSE	ELSA-NSE	6 weeks	370			Fr 28		Fr 25		Fr 25			Fr 22		Fr 20						
17-OH-PROGESTERONE	17-OHP-CT	9 weeks	190				Th 10				Th 24			Th 05					Th 16		
OSTEOCALCIN	OSTEO-RIACT	7 weeks	577		Fr 14			Fr 11		Fr 11		Fr 08			Fr 13						
PROCOLLAGEN III PEPTIDE	OCFK07-PIIIP	6 weeks	300			Fr 28		Fr 25			Fr 25		Fr 22		Fr 20						
PROGESTERONE	PROGEST-RIA	9 weeks	220				Th 10				Th 24			Th 05					Th 16		
PSA	PSA-RIACT	5 weeks	370	We 05			We 02		We 02		We 30		We 27		We 25						
FREE PSA	FPSA-RIACT	5 weeks	185	We 05			We 02		We 02		We 30		We 27		We 25						
PTH	ELSA-PTH	6 weeks	555		We 19		We 16			We 16		We 13		We 11	We 25						
RENIN	RENINE	6 weeks	360		Mo 17			Mo 14			Mo 14		Mo 11		Mo 09	Mo 23					
SHBG	SHBG-RIACT	11 weeks	185				Fr 11					Fr 08			Fr 13						
TBG	OCRU03-TBG	13 weeks	54		We 12					We 09				We 04							
TESTOSTERONE	TESTO-CT2	13 weeks	200				Fr 11					Fr 08			Mo 30						
THYROGLOBULIN	THYRO	6 weeks	481		We 19		We 16			We 16		We 13			We 18						
T3/T4	OCPE07-T3	10 weeks	150		We 12		We 09		We 09			We 06		We 04							
	OCPE57-T3	10 weeks	750		We 12		We 09		We 09			We 06		We 04							
	OCPG07-T4	10 weeks	150		We 12		We 09		We 09			We 06		We 04							
	OCPG57-T4	10 weeks	750		We 12		We 09		We 09			We 06		We 04							
TRYPSIN	OCFE07-TRYPS	8 weeks	90			We 26		We 23		We 23			Th 21		We 18						
TSH	OCPL07-TSH	8 weeks	300	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						
	OCPL57-TSH	8 weeks	1500	Fr 07			Fr 04			Fr 04		Fr 01	Fr 29		Mo 30						

### FOR MORE INFORMATION

■ **Cisbio Bioassays** - Parc Marcel Boiteux, BP 84175, 30200 Codolet - France / **Europe and other countries:** +33 169 857 253 or +33 169 857 546 **Hotline:** +33 466 796 819 **Email:** iva@cisbio.com [www.cisbio.com](http://www.cisbio.com)

# 2022 AVAILABILITY SCHEDULE by product family / grouped availability

In-Vitro RIA Products - See below products concerned by grouped availability.

At least 5 weeks of validity for all products (Except IRMA-HCT (4 weeks)) - An overlap of 9 days between 2 lots.

**BANK HOLIDAYS IN 2022:** 1<sup>st</sup> January, 18<sup>th</sup> April, 1<sup>st</sup> May, 8<sup>th</sup> May, 26<sup>th</sup> May, 6<sup>th</sup> June, 14<sup>th</sup> July, 15<sup>th</sup> August, 1<sup>st</sup> November, 11<sup>th</sup> November, 25<sup>th</sup> December.  
Please order one day before normal ordering day

**ORDERS:** Phone: +33 169 857 253 or +33 169 857 546  
Fax: +33 169 857 521

JANUARY		FEBRUARY		MARCH			APRIL		MAY		JUNE	
05	19	02	16	02	16	30	13	27	11	25	08	22
OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3	OCFH57-FT3
IRMA-CPEP	IRMA-CPEP	CORT-CT2	IRMA-CPEP	IRMA-CPEP	OCPL07-TSH	IRMA-CPEP	IRMA-CPEP	CORT-CT2	IRMA-CPEP	IRMA-CPEP	OCPL07-TSH	CORT-CT2
CORT-CT2	OCPL07-TSH	IRMA-HCT	OCPL07-TSH	CORT-CT2	OCPL57-TSH	CORT-CT2	OCPL07-TSH	IRMA-HCT	OCPL07-TSH	CORT-CT2	OCPL57-TSH	TESTO-CT2
IRMA-HCT	OCPL57-TSH	HGH-RIACT	OCPL57-TSH	IRMA-HCT	CORT-CT2	IRMA-HCT	OCPL57-TSH	HGH-RIACT	OCPL57-TSH	HGH-RIACT	CORT-CT2	SHBG-RIACT
HGH-RIACT	CORT-CT2	IGF1-RIACT	CORT-CT2	HGH-RIACT	IRMA-HCT	HGH-RIACT	CORT-CT2	OCPM03-HCG	CORT-CT2	THYRO	IGF1-RIACT	TGAB I STEP
OCPM03-HCG	IRMA-HCT	BI-INS-IRMA	IRMA-HCT	OCPM03-HCG	HGH-RIACT	IGF1-RIACT	IRMA-HCT	BI-INS-IRMA	IRMA-HCT	OSTEO-RIACT	ALDO-RIACT	REN-CT2P
BI-INS-IRMA	HGH-RIACT	ALDO-RIACT	HGH-RIACT	BI-INS-IRMA	THYRO	BI-INS-IRMA	HGH-RIACT	TESTO-CT2	HGH-RIACT	IGF1-RIACT	TESTO-CT2	PROGEST-RIA
TESTO-CT2	THYRO	TESTO-CT2	THYRO	TESTO-CT2	OSTEO-RIACT	ALDO-RIACT	THYRO	OCPE07-T3	OCPM03-HCG	BI-INS-IRMA	OCPE07-T3	
OCPE07-T3	OSTEO-RIACT	OCPE07-T3	OSTEO-RIACT	OCPE07-T3	OCPM03-HCG	TESTO-CT2	OSTEO-RIACT	OCPE57-T3	IGF1-RIACT	ALDO-RIACT	OCPE57-T3	
OCPE57-T3	OCPM03-HCG	OCPE57-T3	OCPM03-HCG	OCPE57-T3	IGF1-RIACT	OCPE07-T3	OCPM03-HCG	OCPG07-T4	ALDO-RIACT	RENINE	OCPG07-T4	
OCPG07-T4	IGF1-RIACT	OCPG07-T4	IGF1-RIACT	OCPG07-T4	ALDO-RIACT	OCPE57-T3	IGF1-RIACT	OCPG57-T4	RENINE	TESTO-CT2	OCPG57-T4	
OCPG57-T4	ALDO-RIACT	OCPG57-T4	ALDO-RIACT	OCPG57-T4	RENINE	OCPG07-T4	ALDO-RIACT	SHBG-RIACT	TESTO-CT2	OCPE07-T3	SHBG-RIACT	
SHBG-RIACT	RENINE	CGA-RIACT	RENINE	SHBG-RIACT	TESTO-CT2	OCPG57-T4	RENINE	CGA-RIACT	OCPE07-T3	OCPE57-T3	TGAB I STEP	
CGA-RIACT	TESTO-CT2	OCFK07-PIIIP	TESTO-CT2	CGA-RIACT	OCPE07-T3	CGA-RIACT	TESTO-CT2	OCFK07-PIIIP	OCPE57-T3	OCPG07-T4	PROGEST-RIA	
OCFM07-CA50	OCPE07-T3	TGAB I STEP	OCPE07-T3	OCFM07-CA50	OCPE57-T3	OCFM07-CA50	OCPE07-T3	TGAB I STEP	OCPG07-T4	OCPG57-T4		
OCFK07-PIIIP	OCPE57-T3	REN-CT2P	OCPE57-T3	OCFK07-PIIIP	OCPG07-T4	OCFK07-PIIIP	OCPE57-T3	REN-CT2P	OCPG57-T4	SHBG-RIACT		
TGAB I STEP	OCPG07-T4	PROGEST-RIA	OCPG07-T4	TGAB I STEP	OCPG57-T4	TGAB I STEP	OCPG07-T4	PROGEST-RIA	SHBG-RIACT	CGA-RIACT		
REN-CT2P	OCPG57-T4		OCPG57-T4	REN-CT2P	SHBG-RIACT	REN-CT2P	OCPG57-T4		CGA-RIACT	OCFM07-CA50		
PROGEST-RIA	SHBG-RIACT		SHBG-RIACT		CGA-RIACT	PROGEST-RIA	SHBG-RIACT		OCFM07-CA50	OCFK07-PIIIP		
	CGA-RIACT		OCFM07-CA50		TGAB I STEP		OCFM07-CA50		TGAB I STEP	TGAB I STEP		
	OCFM07-CA50		TGAB I STEP		PROGEST-RIA		TGAB I STEP		PROGEST-RIA	REN-CT2P		
	TGAB I STEP		PROGEST-RIA									