

**NEW**  
2021 ANNUAL SCHEDULE

# Annual Availability Schedule 2021

In-Vitro RIA Products

# 2021 - ANNUAL AVAILABILITY SCHEDULE by product

## In-Vitro RIA Products

PARAMETER	KIT REFERENCE	SHELF-LIFE (weeks)	ACT (kBq)	JANUARY			FEBRUARY			MARCH			APRIL			MAY		
				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
ACTH	ELSA-ACTH	7 weeks	592			We 20		We 17			We 17		We 14			We 12		
AFP	ELSA2-AFP	7 weeks	277.5			We 27			We 24		We 24			We 21		We 19		
	OCFA07-AFP	10 weeks	300				We 03				We 31						We 26	
ALDOSTERONE	ALDO-RIACT	9 weeks	111					Mo 15					Mo 12					
ANDROSTENEDIONE	R-GM-100	9 weeks	115		We 13			We 10		We 10		Th 08			We 05			
ANGIOTENSIN I	REN-CT2P	7 weeks	150			Th 28			Th 25		Th 25			Th 22		Th 20		
ANTI-ACETYLCHOLINE RECEPTOR AUTOANTIBODIES	ACHR-AB	6 weeks	40		We 13		We 03			We 03		We 31		We 28			We 26	
ANTI-INSULIN ANTIBODIES	AAI	6 weeks	52	Tu 05			Mo 01			Mo 01		Mo 29		Mo 26			Tu 25	
ANTI-THYROGLOBULINE ANTIBODIES	TGAB I STEP	8 weeks	137		Th 14			Th 11			Th 11		Fr 09			Th 06		
ANTI-TPO AUTOANTIBODIES	TPO-AB-CT	10 weeks	40		We 13		We 03			We 03		We 31		We 28			We 26	
ANTI-TSH RECEPTOR AUTOANTIBODIES	TR-AB-CT	6 weeks	100			We 27			We 24			We 24		We 21		We 19		
ANTI-VGKC AUTOANTIBODIES	VGK/25	8 weeks	30					We 10					We 07					
ANTI-VGCC AUTOANTIBODIES	VGC/25	9 weeks	12					We 10					We 07					
CALCITONIN	IRMA-HCT	4 weeks	611	We 06		We 20	We 03	We 17		We 03	We 17	We 31		We 14	We 28		We 12	We 26
CA 125	ELSA-CA125II	7 weeks	481	We 06			We 03			We 03		We 31		We 28				We 26
CA 15-3	ELSA-CA15-3	7 weeks	555			We 27			We 24			We 24		We 21		We 19		
CA 19-9	ELSA-CA19-9	7 weeks	370			We 27			We 24			We 24		We 21		We 19		
CA 50	OCFM07-CA50	7 weeks	300				We 03				We 17			We 28				
CA72-4	ELSA-CA72-4	7 weeks	370			We 27			We 24			We 24		We 21		We 19		
CEA	CEA-RIACT	6 weeks	350	Tu 05			Mo 01			Mo 01		Mo 29		Mo 26			Tu 25	
CHROMOGRANIN A	CGA-RIACT	9 weeks	407				We 03					We 31					We 26	
CORTISOL	CORT-CT2	9 weeks	250		We 13			We 10			We 10		Th 08			We 05		
C-PEPTIDE	IRMA-CPEP	8 weeks	460				We 03				We 17			We 28				
CYFRA-21-1	ELSA-CYFRA-	7 weeks	277.5	Mo 04		Fr 29			Fr 26			Fr 26		Fr 23			Fr 21	
ESTRADIOL	ESTR-CTRIA	9 weeks	93		Mo 18						Mo 15						Mo 10	
	ESTR-US-CT	7 weeks	85		Fr 15			Fr 12			Fr 12		Fr 09			Fr 07		
FT3/FT4	OCFH07-FT3	9 weeks	150	Fr 08			Fr 05			Fr 05			Fr 02		Fr 30		Fr 28	
	OCFH57-FT3	9 weeks	750	Fr 08			Fr 05			Fr 05			Fr 02		Fr 30		Fr 28	
	OCFD03-FT4	9 weeks	150	Fr 08			Fr 05			Fr 05			Fr 02		Fr 30		Fr 28	
	OCFD03-FT4-K	9 weeks	150	Fr 08			Fr 05			Fr 05			Fr 02		Fr 30		Fr 28	
	OCFD53-FT4	9 weeks	750	Fr 08			Fr 05			Fr 05			Fr 02		Fr 30		Fr 28	
GASTRIN	GASK-PR	7 weeks	37		We 13			We 10			We 10		Th 08			We 05		
HCG - Free $\beta$ chains	ELSA-FBHC	7 weeks	555			We 27			We 24			We 24		We 21		We 19		
HCG	OCPM03-HCG	10 weeks	300			We 20					We 17					We 12		
HGH	HGH-RIACT	7 weeks	308			We 20		We 17			We 17		We 14			We 12		
IGF1	IGF1-RIACT	10 weeks	370					Mo 15					Mo 12					
INSULIN	BI-INS-IRMA	6 weeks	460	Mo 04		Fr 29			Fr 26			Fr 26		Fr 23			Fr 21	
MuSK	MSK/TR	6 weeks	50			We 27			We 24			We 24		We 21		We 19		
NSE	ELSA-NSE	6 weeks	370	Mo 04		Fr 29			Fr 26			Fr 26		Fr 23			Fr 21	
17-OH-PROGESTERONE	17-OHP-CT	9 weeks	190			Th 28					Th 11			Th 22				
OSTEOCALCIN	OSTEO-RIACT	7 weeks	577		Fr 15			Fr 12			Fr 12		Fr 09			Fr 07		
PROCOLLAGEN III PEPTIDE	OCFK07-PIIP	6 weeks	300	Mo 04		Fr 29			Fr 26			Fr 26		Fr 23			Fr 21	
PROGESTERONE	PROGEST-RIA	9 weeks	220			Th 28					Th 11			Th 22				
PSA	PSA-RIACT	5 weeks	370	We 06			We 03			We 03		We 31		We 28			We 26	
FREE PSA	FPSA-RIACT	5 weeks	185	We 06			We 03			We 03		We 31		We 28			We 26	
PTH	ELSA-PTH	6 weeks	555			We 20		We 17			We 17		We 14			We 12		
RENIN	RENINE	6 weeks	360		Mo 18			Mo 15			Mo 15		Mo 12			Mo 10		
SHBG	SHBG-RIACT	11 weeks	185		Fr 15						Fr 12					Fr 07		
TBG	OCRU03-TBG	13 weeks	54					We 10					Th 08					
TESTOSTERONE	TESTO-CT2	13 weeks	200		Fr 15						Fr 12					Fr 07		
THYROGLOBULIN	THYRO	6 weeks	481			We 20		We 17			We 17		We 14			We 12		
T3/T4	OCPE07-T3	10 weeks	150		We 13			We 10			We 10		Th 08			We 05		
	OCPE57-T3	10 weeks	750		We 13			We 10			We 10		Th 08			We 05		
	OCPG07-T4	10 weeks	150		We 13			We 10			We 10		Th 08			We 05		
	OCPG57-T4	10 weeks	750		We 13			We 10			We 10		Th 08			We 05		
TRYPsin	OCFE07-TRYPS	8 weeks	90			We 27			We 24			We 24		We 21		We 19		
TSH	OCPL07-TSH	8 weeks	300	Fr 08			Fr 05			Fr 05			Fr 02	Fr 30			Fr 28	
	OCPL57-TSH	8 weeks	1500	Fr 08			Fr 05			Fr 05			Fr 02	Fr 30			Fr 28	

### 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)

**BANK HOLIDAYS IN 2021:**

1<sup>st</sup> January, 5<sup>th</sup> April, 1<sup>st</sup> May, 8<sup>th</sup> May, 13<sup>th</sup> May, 24<sup>th</sup> May, 1<sup>st</sup> June, 14<sup>th</sup> July,  
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



KIT REFERENCE	JUNE			JULY			AUGUST			SEPTEMBER			OCTOBER			NOVEMBER			DECEMBER		
	1-9	10-19	20-30	1-9	10-19	20-31	1-9	10-19	20-31	1-9	10-19	20-30	1-9	10-19	20-31	1-9	10-19	20-30	1-9	10-19	20-31
ELSA-ACTH	We 09			We 07			We 04			We 01		We 29			We 27			We 24			We 22
ELSA2-AFP		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
OCCA07-AFP						We 21				We 15						We 10					
ALDO-RIACT	Mo 07						Mo 02					Mo 27						Mo 22			
R-GM-100	We 02		We 30			We 28			We 25			We 22			We 20		We 17			We 15	
REN-CT2P		Th 17			Fr 16			Th 12		Th 09			Th 07			Fr 05			Th 02		Th 30
ACHR-AB			We 23			We 21		We 18			We 15			We 13		We 10			We 08		
AAI			Mo 21		Mo 19			Mo 16			Mo 13			Mo 11		Mo 08			Mo 06		
TGAB I STEP	Th 03			Th 01		Th 29			Th 26			Th 23			Th 21		Th 18			Th 16	
TPO-AB-CT			We 23			We 21		We 18			We 15			We 13			We 10		We 08		
TR-AB-CT		We 16			Th 15			We 11		We 08			We 06			We 03			We 01		We 22
VGK/25	We 02					We 28						We 22				We 17					
VGC/25	We 02					We 28						We 22				We 17					
IRMA-HCT	We 09		We 23	We 07		We 21	We 04	We 18		We 01	We 15	We 29		We 13	We 27		We 10	We 24	We 08		We 22
ELSA-CA125II			We 23			We 21		We 18			We 15			We 13			We 10		We 08		
ELSA-CA15-3		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
ELSA-CA19-9		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
OCCM07-CA50	We 09					We 21				We 01				We 13			We 24				
ELSA-CA72-4		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
CEA-RIACT			Mo 21		Mo 19			Mo 16			Mo 13			Mo 11		Mo 08			Mo 06		
CGA-RIACT						We 21					We 15					We 10					
CORT-CT2	We 02		We 30			We 28			We 25			We 22			We 20		We 17			We 15	
IRMA-CPEP	We 09					We 21				We 01				We 13				We 24			
ELSA-CYFRA		Fr 18			Fr 16			Fr 13			Fr 10		Fr 08			Fr 05			Fr 03		Fr 31
ESTR-CTRIA				Mo 05					Mo 30						Mo 25						Mo 20
ESTR-US-CT	Fr 04			Fr 02		Fr 30			Fr 27			Fr 24			Fr 22		Fr 19			Fr 17	
OCCF07-FT3			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
OCCF57-FT3			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
OCCF03-FT4			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
OCCF03-FT4-K			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
OCCF53-FT4			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
GASK-PR	We 02		We 30			We 28			We 25			We 22			We 20		We 17			We 15	
ELSA-FBHC		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
OCCPM03-HCG				We 07						We 01				We 27							We 22
HGH-RIACT	We 09			We 07			We 04			We 01		We 29		We 27			We 24			We 22	
IGF1-RIACT	Mo 07						Mo 02					Mo 27					Mo 22				
BI-INS-IRMA		Fr 18			Fr 16			Fr 13			Fr 10		Fr 08			Fr 05			Fr 03		Fr 31
MSK/TR		We 16			Th 15			We 11		We 08			We 06			We 03			We 01		We 22
ELSA-NSE		Fr 18			Fr 16			Fr 13			Fr 10		Fr 08			Fr 05			Fr 03		Fr 31
17-OHP-CT	Th 03				Fr 16				Th 26				Th 07				Th 18				Th 30
OSTEO-RIACT	Fr 04			Fr 02		Fr 30			Fr 27			Fr 24			Fr 22		Fr 19			Fr 17	
OCCF07-PIIP		Fr 18			Mo 19			Fr 13			Fr 10		Fr 08			Fr 05			Fr 03		Fr 31
PROGEST-RIA	Th 03				Fr 16				Th 26				Th 07				Th 18				Th 30
PSA-RIACT			We 23			We 21		We 18			We 15			We 13			We 10		We 08		
FPSA-RIACT			We 23			We 21		We 18			We 15			We 13			We 10		We 08		
ELSA-PTH	We 09			We 07			We 04			We 01		We 29			We 27		We 24			We 22	
RENINE	Mo 07			Mo 05			Mo 02		Mo 30			Mo 27			Mo 25			Mo 22			Mo 20
SHBG-RIACT				Fr 02					Fr 27					Fr 22						Fr 17	
OCRU03-TBG	We 02					We 28						We 22				We 17					
TESTO-CT2				Fr 02					Fr 27						Fr 22					Fr 17	
THYRO	We 09			We 07			We 04			We 01		We 29		We 27			We 24			We 22	
OCCPE07-T3	We 02		We 30			We 28			We 25			We 22		We 20		We 17			We 15		
OCCPE57-T3	We 02		We 30			We 28			We 25			We 22		We 20		We 17			We 15		
OCCPG07-T4	We 02		We 30			We 28			We 25			We 22		We 20		We 17			We 15		
OCCPG57-T4	We 02		We 30			We 28			We 25			We 22		We 20		We 17			We 15		
OCCFE07-TRYPS		We 16			Th 15			We 11		We 08			We 06			Th 04			We 01		We 29
OCCPL07-TSH			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	
OCCPL57-TSH			Fr 25			Fr 23			Fr 20		Fr 17			Fr 15			Mo 15			Fr 10	





**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](mailto:iva@cisbio.com) **www.cisbio.com**