

**NEW**  
2019 ANNUAL SCHEDULE

# Annual Availability Schedule 2019

In-Vitro RIA Products

# 2019 - 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 23			We 20				We 17			We 15		
AFP	ELSA2-AFP	7 weeks	277.5	Th 03		We 30			We 27		We 27			Th 25			We 22	
	OCFA07-AFP	10 weeks	300				We 06					We 03					We 29	
ALDOSTERONE	ALDO-RIACT	9 weeks	111					Mo 18					Mo 15					
ANDROSTENEDIONE	R-GM-100	9 weeks	115		We 16			We 13		We 13			We 10		Th 09			
ANGIOTENSIN I ANTI-ACETYLCHOLINE	REN-CT2	7 weeks	148	We 09			We 06			We 06		We 03			Th 02		We 29	
RECEPTOR AUTOANTIBODIES	ACHR-AB	6 weeks	40	We 09			Tu 05			Tu 05		Tu 02		Tu 30			Tu 28	
ANTI-INSULIN ANTIBODIES	AAI	6 weeks	52	Mo 07			Mo 04			Mo 04		Mo 01		Mo 29			Mo 27	
ANTI-THYROGLOBULINE ANTIBODIES	TGAB I STEP	7 weeks	137		Th 17			Th 14		Th 14			Th 11			Th 16		
ANTI-TPO AUTOANTIBODIES	TPO-AB-CT	10 weeks	40		Tu 15		Tu 05			Tu 05		Tu 02		Tu 30			Tu 28	
ANTI-TSH RECEPTOR AUTOANTIBODIES	TR-AB-CT	6 weeks	100		Tu 15	Tu 29			Tu 26					We 24			Tu 21	
ANTI-VGKC AUTOANTIBODIES	VGK/25	8 weeks	30					Tu 12				Tu 09						
ANTI-VGCC AUTOANTIBODIES	VGC/25	9 weeks	12					Tu 12				Tu 09						
CALCITONIN	IRMA-HCT	4 weeks	611	We 09		We 23	We 06		We 20	We 06		We 20	We 03	We 17	Th 02	Fr 17	We 29	
CA 125	ELSA-CA125II	7 weeks	481	We 09			We 06			We 06		We 03			Th 02		We 29	
CA 15-3	ELSA-CA15-3	7 weeks	555	Th 03		We 30			We 27			We 27		Th 25			We 22	
CA 19-9	ELSA-CA19-9	7 weeks	370	Th 03		We 30			We 27			We 27		Th 25			We 22	
CA 50	OCFM07-CA50	7 weeks	300	We 09					We 20			We 03				We 15		
CA72-4	ELSA-CA72-4	7 weeks	370	Th 03		We 30			We 27			We 27		Th 25			We 22	
CEA	CEA-RIACT	6 weeks	350	Mo 07			Mo 04			Mo 04		Mo 01		Mo 29			Mo 27	
CHROMOGRANIN A	CGA-RIACT	9 weeks	407				We 06					We 03					We 29	
CORTISOL	CORT-CT2	9 weeks	250		We 16			We 13		We 13			We 10		Th 09			
C-PEPTIDE	IRMA-CPEP	8 weeks	460	We 09					We 20			We 03				We 15		
CYFRA-21-1	ELSA-CYFRA-	7 weeks	277.5	Fr 04			Fr 01			Fr 01		Fr 29		Fr 26			Fr 24	
ESTRADIOL	ESTR-CTRIA	9 weeks	93			Mo 21				Mo 18						Mo 13		
	ESTR-US-CT	7 weeks	85		Fr 18			Fr 15		Fr 15			Fr 12			Mo 13		
FT3/FT4	OCFH07-FT3	9 weeks	150		Fr 11		Fr 08			Fr 08		Fr 05			Mo 06			
	OCFH57-FT3	9 weeks	750		Fr 11		Fr 08			Fr 08		Fr 05			Mo 06			
	OCFD03-FT4	9 weeks	150		Fr 11		Fr 08			Fr 08		Fr 05			Mo 06			
	OCFD03-FT4-K	9 weeks	150		Fr 11		Fr 08			Fr 08		Fr 05			Mo 06			
	OCFD53-FT4	9 weeks	750		Fr 11		Fr 08			Fr 08		Fr 05			Mo 06			
GASTRIN	GASK-PR	7 weeks	37		We 16			We 13		We 13			We 10			Fr 10		
HCG - Free b chains	ELSA-FBHC	7 weeks	555	Th 03		We 30			We 27		We 27		Th 25			We 22		
HCG	OCPM03-HCG	10 weeks	300			We 23					We 20					We 15		
HGH	HGH-RIACT	7 weeks	308			We 23			We 20		We 20		We 17			We 15		
IGF1	IGF1-RIACT	10 weeks	370					Mo 18					Mo 15					
INSULIN	BI-INS-IRMA	6 weeks	460	Fr 04			Fr 01			Fr 01		Fr 29		Fr 26			Fr 24	
MuSK	MSK/TR	6 weeks	50		Tu 15	Tu 29			Tu 26			Tu 26		We 24			Tu 21	
NSE	ELSA-NSE	6 weeks	370	Fr 04			Fr 01			Fr 01		Fr 29		Fr 26			Fr 24	
17-OH-PROGESTERONE	OHP-CT	7 weeks	97		Mo 14			Mo 11		Mo 11		Mo 08			Mo 06			
OSTEOCALCIN	OSTEO-RIACT	7 weeks	577		Fr 18			Fr 15		Fr 15			Fr 12			Mo 13		
PROCOLLAGEN III PEPTIDE	OCFK07-PIIP	6 weeks	300	Fr 04			Fr 01			Fr 01		Fr 29		Fr 26			Fr 24	
PROGESTERONE	PROGEST-RIA	8 weeks	220					Th 14				Th 28				Fr 10		
PSA	PSA-RIACT	5 weeks	370	We 09			We 06			We 06		We 03			Th 02		We 29	
FREE PSA	FPSA-RIACT	5 weeks	185	We 09			We 06			We 06		We 03			Th 02		We 29	
PTH	ELSA-PTH	6 weeks	555			We 23			We 20		We 20		We 17			Fr 17		
RENIN	RENINE	6 weeks	360			Mo 21		Mo 18		Mo 18			Mo 15			Mo 13		
SHBG	SHBG-RIACT	11 weeks	185		Fr 18					Fr 15						Mo 13		
TBG	OCRU03-TBG	13 weeks	54					We 13					We 10					
TESTOSTERONE	TESTO-CT2	13 weeks	200		Fr 18					Fr 15						Fr 10		
THYROGLOBULIN	THYRO	6 weeks	481			We 23			We 20		We 20		We 17			Fr 17		
T3/T4	OCPE07-T3	10 weeks	150		We 16			We 13		We 13		We 10		Th 09				
	OCPE57-T3	10 weeks	750		We 16			We 13		We 13		We 10		Th 09				
	OCPG07-T4	10 weeks	150		We 16			We 13		We 13		We 10		Th 09				
	OCPG57-T4	10 weeks	750		We 16			We 13		We 13		We 10		Th 09				
TRYPsin	OCFE07-TRYPS	8 weeks	90	Th 03		We 30			We 27		We 27		Th 25				We 22	
TSH	OCPL07-TSH	8 weeks	300		Fr 18			Fr 15		Fr 15			Fr 12			Mo 13		
	OCPL57-TSH	8 weeks	1500		Fr 18			Fr 15		Fr 15			Fr 12			Mo 13		

### 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: +33 466 796 787 Email: iva@cisbio.com [www.cisbio.com](http://www.cisbio.com)

**BANK HOLIDAYS IN 2019:**

1<sup>st</sup> January, 22<sup>nd</sup> April, 1<sup>st</sup> May, 8<sup>th</sup> May, 30<sup>th</sup> May, 10<sup>th</sup> June, 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



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 12			We 10		We 07			We 04			We 02		We 30			We 27			Th 26	
ELSA2-AFP		We 19			We 17			We 14			We 11		We 09			We 06				We 04		
OCFA07-AFP						We 24					We 18						Th 14					
ALDO-RIACT		Tu 11					Mo 05					Mo 30						Mo 25				
R-GM-100	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
REN-CT2			We 26			We 24			We 21		We 18			We 16			Th 14				We 11	
ACHR-AB			Tu 25			Tu 23			Tu 20		Tu 17			Tu 15			We 13				Tu 10	
AAI			Mo 24			Mo 22		Mo 19			Mo 16			Mo 14			Tu 12			Mo 09		
TGAB I STEP		Th 13			Th 11			Fr 16			Th 19			Th 17			Fr 15				Th 12	
TPO-AB-CT			Tu 25			Tu 23			Tu 20		Tu 17			Tu 15			We 13				Tu 10	
TR-AB-CT		Tu 18			Tu 16			Tu 13			Tu 10			Tu 08			Tu 05			Tu 03		
VGK/25	Tu 04					Tu 30						Tu 24					Tu 19					
VGC/25	Tu 04					Tu 30						Tu 24					Tu 19					
IRMA-HCT		We 12	We 26		We 10	We 24	We 07		Fr 23	We 04	We 18		We 02	We 16	We 30		Th 14	We 27		We 11	Th 26	
ELSA-CA125II			We 26			We 24			We 21		We 18			We 16			Th 14				We 11	
ELSA-CA15-3		We 19			We 17			We 14			We 11		We 09			We 06				We 04		
ELSA-CA19-9		We 19			We 17			We 14			We 11		We 09			We 06				We 04		
OCFM07-CA50			We 26				We 07				We 18				We 30						We 11	
ELSA-CA72-4		We 19			We 17			We 14			We 11		We 09			We 06				We 04		
CEA-RIACT			Mo 24			Mo 22		Mo 19			Mo 16			Mo 14			Tu 12			Mo 09		
CGA-RIACT						We 24					We 18						Th 14					
CORT-CT2	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
IRMA-CPEP			We 26				We 07				We 18				We 30						We 11	
ELSA-CYFRA-			Fr 21		Fr 19			Mo 19			Fr 13			Fr 11			Fr 08			Fr 06		
ESTR-CTRIA				Mo 08						Mo 02					Mo 28						Mo 23	
ESTR-US-CT	Fr 07			Fr 05			Fr 02		Fr 30			Fr 27			Fr 25			Fr 22			Fr 20	
OCFH07-FT3	Mo 03		Fr 28			Fr 26			Fr 23			Fr 20		Fr 18			Fr 15				Fr 13	
OCFH57-FT3	Mo 03		Fr 28			Fr 26			Fr 23			Fr 20		Fr 18			Fr 15				Fr 13	
OCFD03-FT4	Mo 03		Fr 28			Fr 26			Fr 23			Fr 20		Fr 18			Fr 15				Fr 13	
OCFD03-FT4-K	Mo 03		Fr 28			Fr 26			Fr 23			Fr 20		Fr 18			Fr 15				Fr 13	
OCFD53-FT4	Mo 03		Fr 28			Fr 26			Fr 23			Fr 20		Fr 18			Fr 15				Fr 13	
GASK-PR	We 05			We 03		We 31		We 28			We 25			We 23			We 20			We 18		
ELSA-FBHCG		We 19			We 17			We 14			We 11		We 09			We 06			We 04			
OCPM03-HCG					We 10					We 04				We 30							Th 26	
HGH-RIACT		We 12			We 10		We 07			We 04			We 02		We 30			We 27			Th 26	
IGF1-RIACT	Fr 07						Mo 05					Mo 30						Mo 25				
BI-INS-IRMA			Fr 21		Fr 19			Mo 19			Fr 13			Fr 11			Fr 08			Fr 06		
MSK/TR		Tu 18			Tu 16			Tu 13			Tu 10			Tu 08			Tu 05			Tu 03		
ELSA-NSE			Fr 21		Fr 19			Mo 19			Fr 13			Fr 11			Fr 08			Fr 06		
OHP-CT	Mo 03			Mo 01		Mo 29			Mo 26			Mo 23			Mo 21		Mo 18			Mo 16		
OSTEO-RIACT	Fr 07			Fr 05			Fr 02		Fr 30			Fr 27			Fr 25			Fr 22			Fr 20	
OCFK07-PIIIP			Fr 21		Fr 19			Mo 19			Fr 13			Fr 11			Fr 08			Fr 06		
PROGEST-RIA			Th 20					Th 01			Th 12				Th 24					Th 05		
PSA-RIACT			We 26			We 24			Th 22		We 18			We 16			Th 14			We 11		
FPSA-RIACT			We 26			We 24			Th 22		We 18			We 16			Th 14			We 11		
ELSA-PTH		We 12			We 10		We 07			We 04			We 02		We 30			We 27			Th 26	
RENINE		Tu 11		Mo 08			Mo 05			Mo 02		Mo 30			Mo 28			Mo 25			Mo 23	
SHBG-RIACT				Fr 05					Fr 30						Fr 25						Fr 20	
OCRU03-TBG	We 05					We 31						We 25					We 20					
TESTO-CT2				Fr 05					Fr 30						Fr 25						Fr 20	
THYRO		We 12			We 10		We 07			We 04			We 02		We 30			We 27			Th 26	
OCPE07-T3	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
OCPE57-T3	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
OCPG07-T4	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
OCPG57-T4	We 05			We 03		We 31			We 28			We 25			We 23			We 20			We 18	
OCFE07-TRYPS		We 19			We 17			We 14			We 11		We 09			We 06			We 04			
OCPL07-TSH	Fr 07			Fr 05			Fr 02		Fr 30			Fr 27			Fr 25			Fr 22			Fr 20	
OCPL57-TSH	Fr 07			Fr 05			Fr 02		Fr 30			Fr 27			Fr 25			Fr 22			Fr 20	

