

**NEW**  
2020 ANNUAL SCHEDULE

# Annual Availability Schedule 2020

In-Vitro RIA Products

# 2020 - 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 22		We 19		We 18			Th 16			We 13		
AFP	ELSA2-AFP	7 weeks	277.5	Th 02		We 29		We 26			We 25			We 22			We 20	
	OCFA07-AFP	10 weeks	300	We 08					We 04					We 29				
ALDOSTERONE	ALDO-RIACT	9 weeks	111			Mo 20				Mo 16						Tu 12		
ANDROSTENEDIONE	R-GM-100	9 weeks	115		We 15			We 12		We 11		We 08			We 06			
ANGIOTENSIN I	REN-CT2P	7 weeks	150	To be communicated														
ANTI-ACETYLCHOLINE RECEPTOR AUTOANTIBODIES	ACHR-AB	6 weeks	40		Tu 14		Tu 04			Tu 03		Tu 31			Tu 28		Tu 26	
ANTI-INSULIN ANTIBODIES	AAI	6 weeks	52	Mo 06			Mo 03			Mo 02		Mo 30			Mo 27		Mo 25	
ANTI-THYROGLOBULINE ANTIBODIES	TGAB I STEP	7 weeks	137	To be communicated														
ANTI-TPO AUTOANTIBODIES	TPO-AB-CT	10 weeks	40		Tu 14		Tu 04			Tu 03		Tu 31			Tu 28		Tu 26	
ANTI-TSH RECEPTOR AUTOANTIBODIES	TR-AB-CT	6 weeks	100		Tu 14	Tu 28			Tu 25			Tu 24			Tu 21		Tu 19	
ANTI-VGKC AUTOANTIBODIES	VGK/25	8 weeks	30			Tu 21					Tu 10					Tu 05		
ANTI-VGCC AUTOANTIBODIES	VGC/25	9 weeks	12			Tu 21					Tu 10					Tu 05		
CALCITONIN	IRMA-HCT	4 weeks	611		Fr 10	We 22	We 05	We 19		We 04	We 18		We 01	Th 16	We 29		We 13	Th 28
CA 125	ELSA-CA125II	7 weeks	481	We 08			We 05			We 04			We 01		We 29			We 27
CA 15-3	ELSA-CA15-3	7 weeks	555	Th 02		We 29		We 26				We 25			We 22			We 20
CA 19-9	ELSA-CA19-9	7 weeks	370	Th 02		We 29		We 26				We 25			We 22			We 20
CA 50	OCFM07-CA50	7 weeks	300			We 22				We 04				Th 16				We 27
CA72-4	ELSA-CA72-4	7 weeks	370	Th 02		We 29		We 26				We 25			We 22			We 20
CEA	CEA-RIACT	6 weeks	350	Mo 06			Mo 03			Mo 02		Mo 30			Mo 27			Mo 25
CHROMOGRANIN A	CGA-RIACT	9 weeks	407	We 08						We 04					We 29			
CORTISOL	CORT-CT2	9 weeks	250		We 15			We 12			We 11		We 08			We 06		
C-PEPTIDE	IRMA-CPEP	8 weeks	460			We 22				We 04				Th 16				We 27
CYFRA-21-1	ELSA-CYFRA-	7 weeks	277.5	Mo 06		Fr 31			Fr 28			Fr 27			Fr 24			Mo 25
ESTRADIOL	ESTR-CTRIA	9 weeks	93					Mo 17						Tu 14				
	ESTR-US-CT	7 weeks	85		Fr 17			Fr 14			Fr 13			Fr 10			Mo 11	
FT3/FT4	OCFH07-FT3	9 weeks	150		Fr 10		Fr 07			Fr 06			Fr 03			Mo 04		Fr 29
	OCFH57-FT3	9 weeks	750		Fr 10		Fr 07			Fr 06			Fr 03			Mo 04		Fr 29
	OCFD03-FT4	9 weeks	150		Fr 10		Fr 07			Fr 06			Fr 03			Mo 04		Fr 29
	OCFD03-FT4-K	9 weeks	150		Fr 10		Fr 07			Fr 06			Fr 03			Mo 04		Fr 29
	OCFD53-FT4	9 weeks	750		Fr 10		Fr 07			Fr 06			Fr 03			Mo 04		Fr 29
GASTRIN	GASK-PR	7 weeks	37		We 15			We 12			We 11		We 08			We 06		
HCG - Free $\beta$ chains	ELSA-FBHC	7 weeks	555	Th 02		We 29			We 26			We 25			We 22			We 20
HCG	OCPM03-HCG	10 weeks	300					We 19						Th 16				
HGH	HGH-RIACT	7 weeks	308			We 22		We 19			We 18			Th 16			We 13	
IGF1	IGF1-RIACT	10 weeks	370			Mo 20					Mo 16						Mo 11	
INSULIN	BI-INS-IRMA	6 weeks	460	Mo 06		Fr 31			Fr 28			Fr 27			Fr 24			Mo 25
MuSK	MSK/TR	6 weeks	50		Tu 14	Tu 28			Tu 25			Tu 24			Tu 21		Tu 19	
NSE	ELSA-NSE	6 weeks	370	Mo 06		Fr 31			Fr 28			Fr 27			Fr 24			Mo 25
17-OH-PROGESTERONE	OHP-CT	7 weeks	97		Mo 13			Mo 10		Mo 09			Mo 06			Mo 04		
OSTEOCALCIN	OSTEO-RIACT	7 weeks	577		Fr 17			Fr 14			Fr 13			Fr 10			Mo 11	
PROCOLLAGEN III PEPTIDE	OCFK07-PIIP	6 weeks	300	Mo 06		Fr 31			Fr 28			Fr 27			Fr 24			Mo 25
PROGESTERONE	PROGEST-RIA	8 weeks	220	To be communicated														
PSA	PSA-RIACT	5 weeks	370	Th 09			We 05			We 04			We 01		We 29			We 27
FREE PSA	FPSA-RIACT	5 weeks	185	Th 09			We 05			We 04			We 01		We 29			We 27
PTH	ELSA-PTH	6 weeks	555			We 22		We 19			We 18			Th 16			Th 14	
RENIN	RENINE	6 weeks	360			Mo 20		Mo 17			Mo 16			Tu 14			Tu 12	
SHBG	SHBG-RIACT	11 weeks	185					Fr 14						Fr 10				
TBG	OCRU03-TBG	13 weeks	54		We 15						We 11					We 06		
TESTOSTERONE	TESTO-CT2	13 weeks	200					Fr 14						Fr 10				
THYROGLOBULIN	THYRO	6 weeks	481			We 22		We 19			We 18			Th 16			Fr 15	
T3/T4	OCPE07-T3	10 weeks	150		We 15			We 12			We 11		We 08			We 06		
	OCPE57-T3	10 weeks	750		We 15			We 12			We 11		We 08			We 06		
	OCPG07-T4	10 weeks	150		We 15			We 12			We 11		We 08			We 06		
	OCPG57-T4	10 weeks	750		We 15			We 12			We 11		We 08			We 06		
TRYPSIN	OCFE07-TRYPS	8 weeks	90	Fr 03		We 29		We 26			We 25			We 22				We 20
TSH	OCPL07-TSH	8 weeks	300		Fr 17			Fr 14			Fr 13			Fr 10			Mo 11	
	OCPL57-TSH	8 weeks	1500		Fr 17			Fr 14			Fr 13			Fr 10			Mo 11	

### 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 2020:**

1<sup>st</sup> January, 13<sup>th</sup> April, 1<sup>st</sup> May, 8<sup>th</sup> May, 21<sup>st</sup> May, 1<sup>st</sup> June, 14<sup>th</sup> July, 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 10		We 08			We 05			We 02		We 30			We 28			We 25			We 23
ELSA2-AFP		We 17			Th 16			We 12		We 09				We 07			We 04			We 02	We 30
OCFA07-AFP			We 24					We 19						We 14						We 09	
ALDO-RIACT				Mo 06					Mo 31						Mo 26						Mo 21
R-GM-100	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
REN-CT2P	To be communicated																				
ACHR-AB			Tu 23			Tu 21		Tu 18			Tu 15			Tu 13			Tu 10			Tu 08	
AAI			Mo 22			Mo 20		Mo 17			Mo 14			Mo 12			Mo 09			Mo 07	
TGAB I STEP	To be communicated																				
TPO-AB-CT			Tu 23			Tu 21		Tu 18			Tu 15			Tu 13			Tu 10			Tu 08	
TR-AB-CT		Tu 16			We 15			Tu 11			Tu 08			Tu 06			Tu 03			Tu 01	
VGK/25			Tu 30						Tu 25						Tu 20						Tu 15
VGC/25			Tu 30						Tu 25						Tu 20						Tu 15
IRMA-HCT		We 10	We 24	We 08		We 22	We 05	We 19		We 02	We 16	We 30		We 14	We 28		Th 12	We 25	We 09		We 23
ELSA-CA125II			We 24			We 22		We 19			We 16			We 14			Th 12		We 09		
ELSA-CA15-3		We 17			Th 16			We 12		We 09			We 07				We 04			We 02	We 30
ELSA-CA19-9		We 17			Th 16			We 12		We 09			We 07				We 04			We 02	We 30
OCFM07-CA50				We 08				We 19				We 30					Th 12				We 23
ELSA-CA72-4		We 17			Th 16			We 12		We 09			We 07				We 04			We 02	We 30
CEA-RIACT			Mo 22			Mo 20		Mo 17			Mo 14			Mo 12			Mo 09			Mo 07	
CGA-RIACT			We 24					We 19						We 14						We 09	
CORT-CT2	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
IRMA-CPEP				We 08				We 19				We 30					Th 12				We 23
ELSA-CYFRA		Fr 19			Fr 17			Fr 14			Fr 11		Fr 09				Fr 06			Fr 04	
ESTR-CTRIA	Mo 08						Mo 03					Mo 28							Mo 23		
ESTR-US-CT	Fr 05			Fr 03		Fr 31			Fr 28			Fr 25			Fr 23				Fr 20		Fr 18
OCFH07-FT3			Fr 26			Fr 24			Fr 21		Fr 18			Fr 16			Fr 13				Fr 11
OCFH57-FT3			Fr 26			Fr 24			Fr 21		Fr 18			Fr 16			Fr 13				Fr 11
OCFD03-FT4			Fr 26			Fr 24			Fr 21		Fr 18			Fr 16			Fr 13				Fr 11
OCFD03-FT4-K			Fr 26			Fr 24			Fr 21		Fr 18			Fr 16			Fr 13				Fr 11
OCFD53-FT4			Fr 26			Fr 24			Fr 21		Fr 18			Fr 16			Fr 13				Fr 11
GASK-PR	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
ELSA-FBHCG		We 17			Th 16			We 12		We 09			We 07				We 04			We 02	We 30
OCPM03-HCG		We 10					We 05					We 30							We 25		
HGH-RIACT		We 10		We 08			We 05			We 02		We 30			We 28			We 25			We 23
IGF1-RIACT				Mo 06					Mo 31						Mo 26						Mo 21
BI-INS-IRMA		Fr 19			Fr 17			Fr 14			Fr 11		Fr 09				Fr 06			Fr 04	
MSK/TR		Tu 16			We 15			Tu 11			Tu 08			Tu 06			Tu 03			Tu 01	
ELSA-NSE		Fr 19			Fr 17			Fr 14			Fr 11		Fr 09				Fr 06			Fr 04	
OHP-CT	Tu 02		Mo 29			Mo 27			Mo 24			Mo 21		Mo 19			Mo 16				Mo 14
OSTEO-RIACT	Fr 05			Fr 03		Fr 31			Fr 28			Fr 25			Fr 23			Fr 20			Fr 18
OCFK07-PIIIP		Fr 19				Mo 20		Fr 14			Fr 11		Fr 09				Fr 06			Fr 04	
PROGEST-RIA	To be communicated																				
PSA-RIACT			We 24			We 22		We 19			We 16			We 14			Th 12			We 09	
FPSA-RIACT			We 24			We 22		We 19			We 16			We 14			Th 12			We 09	
ELSA-PTH		We 10		We 08			We 05			We 02		We 30			We 28			We 25			We 23
RENINE	Mo 08			Mo 06			Mo 03		Mo 31			Mo 28			Mo 26			Mo 23			Mo 21
SHBG-RIACT	Fr 05					Fr 31						Fr 25					Fr 20				
OCRU03-TBG				We 01					We 26						We 21						We 16
TESTO-CT2	Fr 05					Fr 31						Fr 25					Fr 20				
THYRO		We 10		We 08			We 05			We 02		We 30			We 28			We 25			We 23
OCPE07-T3	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
OCPE57-T3	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
OCPG07-T4	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
OCPG57-T4	We 03			We 01		We 29			We 26			We 23			We 21		We 18				We 16
OCFE07-TRYPS		We 17			Th 16			We 12		We 09			We 07				We 04			We 02	Th 31
OCPL07-TSH	Fr 05			Fr 03		Fr 31			Fr 28			Fr 25			Fr 23			Fr 20			Fr 18
OCPL57-TSH	Fr 05			Fr 03		Fr 31			Fr 28			Fr 25			Fr 23			Fr 20			Fr 18





**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**