

NOUVEAU
CALENDRIER DE MISE A DISPOSITION 2019



CALENDRIERS DE MISE A DISPOSITION 2019

Trousses pour radioimmunologie

CALENDRIER 2019 par produits

Trousses pour radioimmunologie

Jours fériés en 2019: 1^{er} janvier, 22 avril, 1^{er} mai, 8 mai, 30 mai, 10 juin, 15 août, 1^{er} novembre, 11 novembre, 25 décembre.

PARAMETRE	REFERENCE TROUSSE	VALIDITE (semaines)	ACT (kBq)	JANVIER			FEVRIER			MARS			AVRIL			MAI		
				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 sem.	592			Me 23			Me 20			Me 20		Me 17			Me 15	
AFP	ELSA2-AFP	7 sem.	277.5	Je 03		Me 30			Me 27			Me 27			Je 25		Me 22	
	OCFA07-AFP	10 sem.	300				Me 06						Me 03				Me 29	
ALDOSTERONE	ALDO-RIACT	9 sem.	111					Lu 18						Lu 15				
ANDROSTENEDIONE	R-GM-100	9 sem.	115		Me 16			Me 13			Me 13			Me 10		Je 09		
ANGIOTENSINE I	REN-CT2	7 sem.	148	Me 09			Me 06			Me 06			Me 03			Je 02	Me 29	
ANTICORPS ANTI-INSULINE	AAI	6 sem.	52	Lu 07			Lu 04			Lu 04			Lu 01		Lu 29		Lu 27	
ANTICORPS ANTI RECEPTEUR ACETYLCHOLINE	ACHR-AB	6 sem.	40	Me 09			Ma 05			Ma 05			Ma 02		Ma 30		Ma 28	
ANTICORPS ANTI- RECEPTEUR TSH	TR-AB-CT	6 sem.	100		Ma 15	Ma 29			Ma 26			Ma 26			Me 24		Ma 21	
ANTICORPS ANTI-THYROGLOBULINE	TGAB I STEP	7 sem.	137		Je 17			Je 14			Je 14			Je 11			Je 16	
ANTICORPS ANTI-THYROPEROXYDASE	TPO-AB-CT	10 sem.	40		Ma 15		Ma 05			Ma 05			Ma 02		Ma 30		Ma 28	
ANTICORPS ANTI-VGKC	VGK/25	8 sem.	30					Ma 12					Ma 09					
ANTICORPS ANTI-VGCC	VGC/25	9 sem.	12					Ma 12					Ma 09					
CA 19-9	ELSA-CA19-9	7 sem.	370	Je 03		Me 30			Me 27			Me 27			Je 25		Me 22	
CA 125	ELSA-CA125II	7 sem.	481	Me 09			Me 06			Me 06			Me 03			Je 02	Me 29	
CA 15-3	ELSA-CA15-3	7 sem.	555	Je 03		Me 30			Me 27			Me 27			Je 25		Me 22	
CA 50	OCFM07-CA50	7 sem.	300	Me 09					Me 20				Me 03				Me 15	
CA 72-4	ELSA-CA72-4	7 sem.	370	Je 03		Me 30			Me 27			Me 27			Je 25		Me 22	
CALCITONINE	IRMA-HCT	4 sem.	611	Me 09		Me 23	Me 06		Me 20	Me 06		Me 20	Me 03	Me 17		Je 02	Ve 17	Me 29
CEA	CEA-RIACT	6 sem.	350	Lu 07			Lu 04			Lu 04			Lu 01		Lu 29		Lu 27	
CHROMOGRANINE A	CGA-RIACT	9 sem.	407				Me 06						Me 03				Me 29	
CORTISOL	CORT-CT2	9 sem.	250		Me 16			Me 13			Me 13			Me 10		Je 09		
C-PEPTIDE	IRMA-CPEP	8 sem.	460	Me 09					Me 20				Me 03				Me 15	
CYFRA-21-1	ELSA-CYFRA-	7 sem.	277.5	Ve 04			Ve 01			Ve 01		Ve 29			Ve 26		Ve 24	
ESTRADIOL	ESTR-CTRIA	9 sem.	93			Lu 21					Lu 18						Lu 13	
	ESTR-US-CT	7 sem.	85		Ve 18			Ve 15			Ve 15			Ve 12			Lu 13	
GASTRINE	GASK-PR	7 sem.	37		Me 16			Me 13			Me 13		Me 10				Ve 10	
HCG (chaines B libres)	ELSA-FBHCg	7 sem.	555	Je 03		Me 30			Me 27			Me 27			Je 25		Me 22	
HCG	OCPM03-HCG	10 sem.	300			Me 23						Me 20					Me 15	
HGH	HGH-RIACT	7 sem.	308			Me 23			Me 20			Me 20		Me 17			Me 15	
IGF1	IGF1-RIACT	10 sem.	370					Lu 18						Lu 15				
INSULIN	BI-INS-IRMA	6 sem.	460	Ve 04			Ve 01			Ve 01		Ve 29			Ve 26		Ve 24	
MuSK	MSK/TR	6 sem.	50		Ma 15	Ma 29			Ma 26			Ma 26			Me 24		Ma 21	
NSE	ELSA-NSE	6 sem.	370	Ve 04			Ve 01			Ve 01		Ve 29			Ve 26		Ve 24	
17-OH-PROGESTERONE	OHP-CT	7 sem.	97		Lu 14			Lu 11			Lu 11		Lu 08			Lu 06		
OSTEOCALCINE	OSTEO-RIACT	7 sem.	577		Ve 18			Ve 15			Ve 15			Ve 12			Lu 13	
PARATHORMONE	ELSA-PTH	6 sem.	555			Me 23			Me 20			Me 20		Me 17			Ve 17	
PEPTIDE PROCOLLAGENE III	OCFK07-PIIIP	6 sem.	300	Ve 04			Ve 01			Ve 01		Ve 29			Ve 26		Ve 24	
PROGESTERONE	PROGEST-RIA	8 sem.	220					Je 14				Je 28					Ve 10	
PSA	PSA-RIACT	5 sem.	370	Me 09			Me 06			Me 06			Me 03			Je 02	Me 29	
PSA LIBRE	FPSA-RIACT	5 sem.	185	Me 09			Me 06			Me 06			Me 03			Je 02	Me 29	
RENINE	RENINE	6 sem.	360			Lu 21		Lu 18			Lu 18			Lu 15			Lu 13	
SHBG	SHBG-RIACT	11 sem.	185		Ve 18						Ve 15						Lu 13	
TBG	OCRU03-TBG	13 sem.	54					Me 13						Me 10				
TESTOSTERONE	TESTO-CT2	13 sem.	200		Ve 18						Ve 15						Ve 10	
THYROGLOBULINE	THYRO	6 sem.	481			Me 23			Me 20			Me 20		Me 17			Ve 17	
THYROXINE	OCPG07-T4	10 sem.	150		Me 16			Me 13			Me 13			Me 10		Je 09		
	OCFD03-FT4	9 sem.	150		Ve 11		Ve 08			Ve 08			Ve 05			Lu 06		
TRIIODOTHYRONINE	OCPE07-T3	10 sem.	150		Me 16			Me 13			Me 13			Me 10		Je 09		
	OCFH07-FT3	9 sem.	150		Ve 11		Ve 08			Ve 08			Ve 05			Lu 06		
TRYPISINE	OCFE07-TRYPS	8 sem.	90	Je 03		Me 30			Me 27			Me 27		Je 25			Me 22	
TSH	OCPL07-TSH	8 sem.	300		Ve 18			Ve 15			Ve 15			Ve 12			Lu 13	

POUR PLUS D'INFORMATIONS

Cisbio Bioassays - Parc Marcel Boiteux, BP 84175, 30200 Codolet - Téléphone : 01 69 85 73 35 Hotline : 04 66 79 67 87 Email : iva@cisbio.com www.cisbio.com

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

