

NOUVEAU
CALENDRIER DE MISE A DISPOSITION 2021



CALENDRIERS DE MISE A DISPOSITION 2021

Trousses pour dosage radioimmunologique

CALENDRIER 2021 par produits

Trousses pour dosage radioimmunologique

Jours fériés en 2021: 1^{er} janvier, 5 avril, 1^{er} mai, 8 mai, 13 mai, 24 mai, 1^{er} juin, 14 juillet, 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 20		Me 17			Me 17		Me 14			Me 12		
AFP	ELSA2-AFP	7 sem.	277.5			Me 27			Me 24			Me 24		Me 21		Me 19		
	OCFA07-AFP	10 sem.	300				Me 03					Me 31					Me 26	
ALDOSTERONE	ALDO-RIACT	9 sem.	111					Lu 15					Lu 12					
ANDROSTENEDIONE	R-GM-100	9 sem.	115		Me 13			Me 10			Me 10		Je 08			Me 05		
ANGIOTENSINE I	REN-CT2P	7 sem.	150			Je 28			Je 25			Je 25		Je 22		Je 20		
ANTICORPS ANTI-INSULINE	AAI	6 sem.	52	Ma 05			Lu 01			Lu 01		Lu 29		Lu 26			Ma 25	
ANTICORPS ANTI RECEPTEUR ACETYLCHOLINE	ACHR-AB	6 sem.	40		Me 13		Me 03			Me 03		Me 31		Me 28			Me 26	
ANTICORPS ANTI- RECEPTEUR TSH	TR-AB-CT	6 sem.	100			Me 27			Me 24			Me 24		Me 21		Me 19		
ANTICORPS ANTI-THYROGLOBULINE	TGAB I STEP	8 sem.	137		Je 14			Je 11			Je 11		Ve 09		Je 06			
ANTICORPS ANTI-THYROPEROXYDASE	TPO-AB-CT	10 sem.	40		Me 13		Me 03			Me 03		Me 31		Me 28			Me 26	
ANTICORPS ANTI-VGKC	VGK/25	8 sem.	30					Me 10					Me 07					
ANTICORPS ANTI-VGCC	VGC/25	9 sem.	12					Me 10					Me 07					
CA 19-9	ELSA-CA19-9	7 sem.	370			Me 27			Me 24			Me 24		Me 21		Me 19		
CA 125	ELSA-CA125II	7 sem.	481	Me 06			Me 03			Me 03		Me 31		Me 28			Me 26	
CA 15-3	ELSA-CA15-3	7 sem.	555			Me 27			Me 24			Me 24		Me 21		Me 19		
CA 50	OCFM07-CA50	7 sem.	300				Me 03				Me 17			Me 28				
CA 72-4	ELSA-CA72-4	7 sem.	370			Me 27			Me 24			Me 24		Me 21		Me 19		
CALCITONINE	IRMA-HCT	4 sem.	611	Me 06		Me 20	Me 03	Me 17		Me 03	Me 17	Me 31		Me 14	Me 28	Me 12	Me 26	
CEA	CEA-RIACT	6 sem.	350	Ma 05			Lu 01			Lu 01		Lu 29		Lu 26			Ma 25	
CHROMOGRANINE A	CGA-RIACT	9 sem.	407				Me 03					Me 31					Me 26	
CORTISOL	CORT-CT2	9 sem.	250		Me 13			Me 10			Me 10		Je 08			Me 05		
C-PEPTIDE	IRMA-CPEP	8 sem.	460				Me 03				Me 17			Me 28				
CYFRA-21-1	ELSA-CYFRA-	7 sem.	277.5	Lu 04		Ve 29			Ve 26			Ve 26		Ve 23			Ve 21	
ESTRADIOL	ESTR-CTRIA	9 sem.	93		Lu 18						Lu 15					Lu 10		
	ESTR-US-CT	7 sem.	85		Ve 15			Ve 12			Ve 12		Ve 09		Ve 07			
GASTRINE	GASK-PR	7 sem.	37		Me 13			Me 10			Me 10		Je 08		Me 05			
HCG (chaines β libres)	ELSA-FBHCG	7 sem.	555			Me 27			Me 24			Me 24		Me 21		Me 19		
HCG	OCPM03-HCG	10 sem.	300			Me 20					Me 17					Me 12		
HGH	HGH-RIACT	7 sem.	308			Me 20		Me 17			Me 17		Me 14			Me 12		
IGF1	IGF1-RIACT	10 sem.	370					Lu 15					Lu 12					
INSULINE	BI-INS-IRMA	6 sem.	460	Lu 04		Ve 29			Ve 26			Ve 26		Ve 23			Ve 21	
MuSK	MSK/TR	6 sem.	50			Me 27			Me 24			Me 24		Me 21		Me 19		
NSE	ELSA-NSE	6 sem.	370	Lu 04		Ve 29			Ve 26			Ve 26		Ve 23			Ve 21	
17-OH-PROGESTERONE	17-OHP-CT	9 sem.	190			Je 28					Je 11			Je 22				
OSTEOCALCINE	OSTEO-RIACT	7 sem.	577		Ve 15			Ve 12			Ve 12		Ve 09		Ve 07			
PARATHORMONE	ELSA-PTH	6 sem.	555			Me 20		Me 17			Me 17		Me 14			Me 12		
PEPTIDE PROCOLLAGENE III	OCFK07-PIIIP	6 sem.	300	Lu 04		Ve 29			Ve 26			Ve 26		Ve 23			Ve 21	
PROGESTERONE	PROGEST-RIA	9 sem.	220			Je 28					Je 11			Je 22				
PSA	PSA-RIACT	5 sem.	370	Me 06			Me 03			Me 03		Me 31		Me 28			Me 26	
PSA LIBRE	FPSA-RIACT	5 sem.	185	Me 06			Me 03			Me 03		Me 31		Me 28			Me 26	
RENINE	RENINE	6 sem.	360		Lu 18			Lu 15			Lu 15		Lu 12			Lu 10		
SHBG	SHBG-RIACT	11 sem.	185		Ve 15						Ve 12				Ve 07			
TBG	OCRU03-TBG	13 sem.	54					Me 10					Je 08					
TESTOSTERONE	TESTO-CT2	13 sem.	200		Ve 15						Ve 12				Ve 07			
THYROGLOBULINE	THYRO	6 sem.	481			Me 20		Me 17			Me 17		Me 14			Me 12		
THYROXINE	OCPG07-T4	10 sem.	150		Me 13			Me 10			Me 10		Je 08		Me 05			
	OCFD03-FT4	9 sem.	150	Ve 08			Ve 05			Ve 05			Ve 02	Ve 30			Ve 28	
TRIIODOTHYRONINE	OCPE07-T3	10 sem.	150		Me 13			Me 10			Me 10		Je 08		Me 05			
	OCFH07-FT3	9 sem.	150	Ve 08			Ve 05			Ve 05			Ve 02	Ve 30			Ve 28	
TRYPISINE	OCFE07-TRYPS	8 sem.	90			Me 27			Me 24			Me 24		Me 21		Me 19		
TSH	OCPL07-TSH	8 sem.	300	Ve 08			Ve 05			Ve 05			Ve 02	Ve 30			Ve 28	

POUR PLUS D'INFORMATIONS

Cisbio Bioassays - Parc Marcel Boiteux, BP 84175, 30200 Codolet FRANCE - Téléphone : 01 69 85 73 35 Hotline : 04 66 79 68 19 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 09			Me 07			Me 04			Me 01		Me 29			Me 27			Me 24			Me 22	
ELSA2-AFP		Me 16			Je 15			Me 11		Me 08				Me 06			Je 04			Me 01		Me 29
OCFA07-AFP						Me 21					Me 15						Me 10					
ALDO-RIACT	Lu 07						Lu 02					Lu 27						Lu 22				
R-GM-100	Me 02		Me 30			Me 28			Me 25			Me 22			Me 20		Me 17			Me 15		
REN-CT2P		Je 17			Ve 16			Je 12		Je 09				Je 07			Ve 05			Je 02		Je 30
AAI			Lu 21		Lu 19			Lu 16			Lu 13			Lu 11		Lu 08			Lu 06			
ACHR-AB			Me 23			Me 21		Me 18			Me 15			Me 13			Me 10		Me 08			
TR-AB-CT		Me 16			Je 15			Me 11		Me 08			Me 06			Me 03			Me 01			Me 22
TGAB I STEP	Je 03			Je 01		Je 29			Je 26			Je 23			Je 21		Je 18			Je 16		
TPO-AB-CT			Me 23			Me 21		Me 18			Me 15			Me 13			Me 10		Me 08			
VGK/25	Me 02					Me 28						Me 22					Me 17					
VGK/25	Me 02					Me 28						Me 22					Me 17					
ELSA-CA19-9		Me 16			Je 15			Me 11		Me 08			Me 06			Je 04			Me 01			Me 29
ELSA-CA125II			Me 23			Me 21		Me 18			Me 15			Me 13			Me 10		Me 08			
ELSA-CA15-3		Me 16			Je 15			Me 11		Me 08			Me 06			Je 04			Me 01			Me 29
OCFM07-CA50	Me 09					Me 21				Me 01				Me 13				Me 24				
ELSA-CA72-4		Me 16			Je 15			Me 11		Me 08			Me 06			Je 04			Me 01			Me 29
IRMA-HCT	Me 09		Me 23	Me 07		Me 21	Me 04	Me 18		Me 01	Me 15	Me 29		Me 13	Me 27		Me 10	Me 24	Me 08		Me 22	
CEA-RIACT			Lu 21		Lu 19			Lu 16			Lu 13			Lu 11		Lu 08			Lu 06			
CGA-RIACT						Me 21					Me 15						Me 10					
CORT-CT2	Me 02		Me 30			Me 28			Me 25			Me 22			Me 20		Me 17			Me 15		
IRMA-CPEP	Me 09					Me 21				Me 01				Me 13				Me 24				
ELSA-CYFRA-		Ve 18			Ve 16			Ve 13			Ve 10		Ve 08			Ve 05			Ve 03			Ve 31
ESTR-CTRIA				Lu 05					Lu 30						Lu 25							Lu 20
ESTR-US-CT	Ve 04			Ve 02		Ve 30			Ve 27			Ve 24			Ve 22		Ve 19			Ve 17		
GASK-PR	Me 02		Me 30			Me 28			Me 25			Me 22			Me 20		Me 17			Me 15		
ELSA-FBHC		Me 16			Je 15			Me 11		Me 08			Me 06			Je 04			Me 01			Me 29
OCPM03-HCG				Me 07						Me 01					Me 27							Me 22
HGH-RIACT	Me 09			Me 07			Me 04			Me 01		Me 29			Me 27			Me 24				Me 22
IGF1-RIACT	Lu 07						Lu 02					Lu 27						Lu 22				
BI-INS-IRMA		Ve 18			Ve 16			Ve 13			Ve 10		Ve 08			Ve 05			Ve 03			Ve 31
MSK/TR		Me 16			Je 15			Me 11		Me 08			Me 06			Me 03			Me 01			Me 22
ELSA-NSE		Ve 18			Ve 16			Ve 13			Ve 10		Ve 08			Ve 05			Ve 03			Ve 31
17-OHP-CT	Je 03				Ve 16				Je 26				Je 07			Je 18						Je 30
OSTEO-RIACT	Ve 04			Ve 02		Ve 30			Ve 27			Ve 24			Ve 22		Ve 19			Ve 17		
ELSA-PTH	Me 09			Me 07			Me 04			Me 01		Me 29			Me 27			Me 24				Me 22
OCFK07-PIIIP		Ve 18			Lu 19			Ve 13			Ve 10		Ve 08			Ve 05			Ve 03			Ve 31
PROGEST-RIA	Je 03				Ve 16				Je 26				Je 07			Je 18						Je 30
PSA-RIACT			Me 23			Me 21		Me 18			Me 15			Me 13			Me 10		Me 08			
FPSA-RIACT			Me 23			Me 21		Me 18			Me 15			Me 13			Me 10		Me 08			
RENINE	Lu 07			Lu 05			Lu 02		Lu 30			Lu 27			Lu 25			Lu 22				Lu 20
SHBG-RIACT				Ve 02					Ve 27					Ve 22						Ve 17		
OCRU03-TBG	Me 02					Me 28						Me 22				Me 17						
TESTO-CT2				Ve 02					Ve 27						Ve 22					Ve 17		
THYRO	Me 09			Me 07			Me 04			Me 01		Me 29			Me 27			Me 24				Me 22
OCPG07-T4	Me 02		Me 30			Me 28			Me 25			Me 22			Me 20		Me 17			Me 15		
OCFD03-FT4			Ve 25			Ve 23			Ve 20		Ve 17			Ve 15			Lu 15			Ve 10		
OCPE07-T3	Me 02		Me 30			Me 28		Me 25			Me 22			Me 20		Me 17			Me 15			
OCFH07-FT3			Ve 25			Ve 23			Ve 20		Ve 17			Ve 15			Lu 15			Ve 10		
OCFE07-TRYPS		Me 16			Je 15			Me 11		Me 08			Me 06			Je 04			Me 01			Me 29
OCPL07-TSH			Ve 25			Ve 23			Ve 20		Ve 17			Ve 15			Lu 15			Ve 10		



POUR PLUS D'INFORMATIONS

Cisbio Bioassays - Parc Marcel Boiteux, BP 84175, 30200 Codolet - France

Téléphone : 01 69 85 73 35 Hotline : 04 66 79 68 19 Email : iva@cisbio.com www.cisbio.com