

NOUVEAU
CALENDRIER DE MISE A DISPOSITION 2020



CALENDRIERS DE MISE A DISPOSITION 2020

Trousses pour dosage radioimmunologique

CALENDRIER 2020 par produits

Trousses pour dosage radioimmunologique

Jours fériés en 2020: 1^{er} janvier, 13 avril, 1^{er} mai, 8 mai, 21 mai, 1^{er} juin, 14 juillet, 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 22		Me 19			Me 18		Je 16			Me 13		
AFP	ELSA2-AFP	7 sem.	277.5	Je 02		Me 29			Me 26			Me 25		Me 22			Me 20	
	OCFA07-AFP	10 sem.	300	Me 08						Me 04			Me 29					
ALDOSTERONE	ALDO-RIACT	9 sem.	111			Lu 20					Lu 16					Ma 12		
ANDROSTENEDIONE	R-GM-100	9 sem.	115		Me 15			Me 12			Me 11		Me 08			Me 06		
ANGIOTENSINE I	REN-CT2P	7 sem.	150	A venir														
ANTICORPS ANTI-INSULINE	AAI	6 sem.	52	Lu 06			Lu 03			Lu 02		Lu 30			Lu 27		Lu 25	
ANTICORPS ANTI RECEPTEUR ACETYLCHOLINE	ACHR-AB	6 sem.	40		Ma 14		Ma 04			Ma 03		Ma 31		Ma 28			Ma 26	
ANTICORPS ANTI- RECEPTEUR TSH	TR-AB-CT	6 sem.	100		Ma 14	Ma 28			Ma 25			Ma 24		Ma 21		Ma 19		
ANTICORPS ANTI-THYROGLOBULINE	TGAB I STEP	7 sem.	137	A venir														
ANTICORPS ANTI-THYROPEROXYDASE	TPO-AB-CT	10 sem.	40		Ma 14		Ma 04			Ma 03		Ma 31		Ma 28			Ma 26	
ANTICORPS ANTI-VGKC	VGK/25	8 sem.	30			Ma 21					Ma 10					Ma 05		
ANTICORPS ANTI-VGCC	VGC/25	9 sem.	12			Ma 21					Ma 10					Ma 05		
CA 19-9	ELSA-CA19-9	7 sem.	370	Je 02		Me 29			Me 26			Me 25		Me 22			Me 20	
CA 125	ELSA-CA125II	7 sem.	481	Me 08			Me 05			Me 04			Me 01	Me 29			Me 27	
CA 15-3	ELSA-CA15-3	7 sem.	555	Je 02		Me 29			Me 26			Me 25		Me 22			Me 20	
CA 50	OCFM07-CA50	7 sem.	300			Me 22				Me 04			Je 16				Me 27	
CA 72-4	ELSA-CA72-4	7 sem.	370	Je 02		Me 29			Me 26			Me 25		Me 22			Me 20	
CALCITONINE	IRMA-HCT	4 sem.	611		Ve 10	Me 22	Me 05	Me 19		Me 04	Me 18		Me 01	Je 16	Me 29	Me 13	Je 28	
CEA	CEA-RIACT	6 sem.	350	Lu 06			Lu 03			Lu 02		Lu 30		Lu 27			Lu 25	
CHROMOGRANINE A	CGA-RIACT	9 sem.	407	Me 08						Me 04				Me 29				
CORTISOL	CORT-CT2	9 sem.	250		Me 15			Me 12			Me 11		Me 08			Me 06		
C-PEPTIDE	IRMA-CPEP	8 sem.	460			Me 22				Me 04			Je 16				Me 27	
CYFRA-21-1	ELSA-CYFRA-	7 sem.	277.5	Lu 06		Ve 31			Ve 28			Ve 27		Ve 24			Lu 25	
ESTRADIOL	ESTR-CTRIA	9 sem.	93				Lu 17						Ma 14					
	ESTR-US-CT	7 sem.	85		Ve 17			Ve 14			Ve 13		Ve 10			Lu 11		
GASTRINE	GASK-PR	7 sem.	37		Me 15			Me 12			Me 11		Me 08			Me 06		
HCG (chaines β libres)	ELSA-FBHCG	7 sem.	555	Je 02		Me 29			Me 26			Me 25		Me 22			Me 20	
HCG	OCPM03-HCG	10 sem.	300					Me 19					Je 16					
HGH	HGH-RIACT	7 sem.	308			Me 22		Me 19			Me 18		Je 16			Me 13		
IGF1	IGF1-RIACT	10 sem.	370			Lu 20					Lu 16					Lu 11		
INSULINE	BI-INS-IRMA	6 sem.	460	Lu 06		Ve 31			Ve 28			Ve 27		Ve 24			Lu 25	
MuSK	MSK/TR	6 sem.	50		Ma 14	Ma 28			Ma 25			Ma 24		Ma 21		Ma 19		
NSE	ELSA-NSE	6 sem.	370	Lu 06		Ve 31			Ve 28			Ve 27		Ve 24			Lu 25	
17-OH-PROGESTERONE	OHP-CT	7 sem.	97		Lu 13			Lu 10		Lu 09			Lu 06			Lu 04		
OSTEOCALCINE	OSTEO-RIACT	7 sem.	577		Ve 17			Ve 14			Ve 13		Ve 10			Lu 11		
PARATHORMONE	ELSA-PTH	6 sem.	555			Me 22		Me 19			Me 18		Je 16			Je 14		
PEPTIDE PROCOLLAGENE III	OCFK07-PIIIP	6 sem.	300	Lu 06		Ve 31			Ve 28			Ve 27		Ve 24			Lu 25	
PROGESTERONE	PROGEST-RIA	8 sem.	220	A venir														
PSA	PSA-RIACT	5 sem.	370	Je 09			Me 05			Me 04			Me 01		Me 29			Me 27
PSA LIBRE	FPSA-RIACT	5 sem.	185	Je 09			Me 05			Me 04			Me 01		Me 29			Me 27
RENINE	RENINE	6 sem.	360			Lu 20		Lu 17			Lu 16		Ma 14			Ma 12		
SHBG	SHBG-RIACT	11 sem.	185					Ve 14					Ve 10					
TBG	OCRU03-TBG	13 sem.	54		Me 15						Me 11					Me 06		
TESTOSTERONE	TESTO-CT2	13 sem.	200					Ve 14					Ve 10					
THYROGLOBULINE	THYRO	6 sem.	481			Me 22		Me 19			Me 18		Je 16			Ve 15		
THYROXINE	OCPG07-T4	10 sem.	150		Me 15			Me 12			Me 11		Me 08			Me 06		
	OCFD03-FT4	9 sem.	150		Ve 10		Ve 07			Ve 06			Ve 03		Lu 04		Ve 29	
TRIIODOTHYRONINE	OCPE07-T3	10 sem.	150		Me 15			Me 12			Me 11		Me 08			Me 06		
	OCFH07-FT3	9 sem.	150		Ve 10		Ve 07			Ve 06			Ve 03		Lu 04		Ve 29	
TRYPSINE	OCFE07-TRYPS	8 sem.	90	Ve 03		Me 29			Me 26			Me 25		Me 22			Me 20	
TSH	OCPL07-TSH	8 sem.	300		Ve 17			Ve 14			Ve 13		Ve 10			Lu 11		

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 10		Me 08			Me 05			Me 02		Me 30			Me 28			Me 25			Me 23
ELSA2-AFP		Me 17			Je 16			Me 12		Me 09				Me 07			Me 04			Me 02	Me 30
OCFA07-AFP			Me 24					Me 19						Me 14					Me 09		
ALDO-RIACT				Lu 06					Lu 31						Lu 26						Lu 21
R-GM-100	Me 03			Me 01		Me 29			Me 26			Me 23			Me 21		Me 18			Me 16	
REN-CT2P	A venir																				
AAI			Lu 22			Lu 20		Lu 17			Lu 14			Lu 12		Lu 09			Lu 07		
ACHR-AB			Ma 23			Ma 21		Ma 18			Ma 15			Ma 13			Ma 10		Ma 08		
TR-AB-CT		Ma 16			Me 15			Ma 11		Ma 08			Ma 06			Ma 03			Ma 01		
TGAB I STEP	A venir																				
TPO-AB-CT			Ma 23			Ma 21		Ma 18			Ma 15			Ma 13			Ma 10		Ma 08		
VGK/25			Ma 30						Ma 25					Ma 20						Ma 15	
VGC/25			Ma 30						Ma 25					Ma 20						Ma 15	
ELSA-CA19-9		Me 17			Je 16			Me 12		Me 09			Me 07			Me 04			Me 02		Me 30
ELSA-CA125II			Me 24			Me 22		Me 19			Me 16		Me 14				Je 12		Me 09		
ELSA-CA15-3		Me 17			Je 16			Me 12		Me 09			Me 07			Me 04			Me 02		Me 30
OCFM07-CA50				Me 08				Me 19				Me 30					Je 12				Me 23
ELSA-CA72-4		Me 17			Je 16			Me 12		Me 09			Me 07			Me 04			Me 02		Me 30
IRMA-HCT		Me 10	Me 24	Me 08		Me 22	Me 05	Me 19		Me 02	Me 16	Me 30		Me 14	Me 28		Je 12	Me 25	Me 09		Me 23
CEA-RIACT			Lu 22			Lu 20		Lu 17			Lu 14			Lu 12		Lu 09			Lu 07		
CGA-RIACT			Me 24					Me 19						Me 14					Me 09		
CORT-CT2	Me 03			Me 01		Me 29			Me 26			Me 23			Me 21		Me 18			Me 16	
IRMA-CPEP				Me 08				Me 19				Me 30					Je 12				Me 23
ELSA-CYFRA-		Ve 19			Ve 17			Ve 14			Ve 11		Ve 09			Ve 06			Ve 04		
ESTR-CTRIA	Lu 08						Lu 03					Lu 28							Lu 23		
ESTR-US-CT	Ve 05			Ve 03		Ve 31			Ve 28			Ve 25		Ve 23				Ve 20		Ve 18	
GASK-PR	Me 03			Me 01		Me 29			Me 26			Me 23		Me 21		Me 18				Me 16	
ELSA-FBHC		Me 17			Je 16			Me 12		Me 09			Me 07			Me 04			Me 02		Me 30
OCPM03-HCG		Me 10					Me 05					Me 30						Me 25			
HGH-RIACT		Me 10		Me 08			Me 05			Me 02		Me 30		Me 28				Me 25			Me 23
IGF1-RIACT				Lu 06					Lu 31						Lu 26						Lu 21
BI-INS-IRMA		Ve 19			Ve 17			Ve 14			Ve 11		Ve 09			Ve 06			Ve 04		
MSK/TR		Ma 16			Me 15			Ma 11		Ma 08			Ma 06			Ma 03			Ma 01		
ELSA-NSE		Ve 19			Ve 17			Ve 14			Ve 11		Ve 09			Ve 06			Ve 04		
OHP-CT	Ma 02		Lu 29			Lu 27			Lu 24			Lu 21		Lu 19			Lu 16			Lu 14	
OSTEO-RIACT	Ve 05			Ve 03		Ve 31			Ve 28			Ve 25		Ve 23				Ve 20		Ve 18	
ELSA-PTH		Me 10		Me 08			Me 05			Me 02		Me 30		Me 28				Me 25			Me 23
OCFK07-PIIIP		Ve 19				Lu 20		Ve 14			Ve 11		Ve 09			Ve 06			Ve 04		
PROGEST-RIA	A venir																				
PSA-RIACT			Me 24			Me 22		Me 19			Me 16			Me 14			Je 12		Me 09		
FPSA-RIACT			Me 24			Me 22		Me 19			Me 16			Me 14			Je 12		Me 09		
RENINE	Lu 08			Lu 06			Lu 03		Lu 31			Lu 28		Lu 26					Lu 23		Lu 21
SHBG-RIACT	Ve 05					Ve 31						Ve 25						Ve 20			
OCRU03-TBG				Me 01					Me 26					Me 21						Me 16	
TESTO-CT2	Ve 05					Ve 31						Ve 25						Ve 20			
THYRO		Me 10		Me 08			Me 05			Me 02		Me 30		Me 28				Me 25			Me 23
OCPG07-T4	Me 03			Me 01		Me 29			Me 26			Me 23		Me 21		Me 18				Me 16	
OCFD03-FT4			Ve 26			Ve 24			Ve 21		Ve 18			Ve 16			Ve 13			Ve 11	
OCPE07-T3	Me 03			Me 01		Me 29			Me 26			Me 23		Me 21		Me 18				Me 16	
OCFH07-FT3			Ve 26			Ve 24			Ve 21		Ve 18			Ve 16			Ve 13			Ve 11	
OCFE07-TRYPS		Me 17			Je 16			Me 12		Me 09			Me 07			Me 04			Me 02		Je 31
OCPL07-TSH	Ve 05			Ve 03		Ve 31			Ve 28			Ve 25		Ve 23			Ve 20			Ve 18	

CALENDRIER 2020 par famille de produits / disponibilités groupées



Trousses pour dosage radioimmunologique - Ci-dessous les produits concernés par les disponibilités groupées.

Au moins 5 semaines de validité pour tous les produits (Excepté IRMA-HCT (4 semaines)) - Un recouvrement de 9 jours entre chaque lot.

Jours fériés en 2020: 1^{er} janvier, 13 avril, 1^{er} mai, 8 mai, 21 mai, 1^{er} juin, 14 juillet, 11 novembre, 25 décembre.

Service commercial / Service COMMANDE: Téléphone : 01 69 85 73 35
Fax : 01 69 85 73 65

JANVIER		FEVRIER		MARS		AVRIL			MAI		JUIN		JUILLET		AOÛT		SEPTEMBRE			OCTOBRE		NOVEMBRE		DECEMBRE	
08	22	05	19	04	18	01	15	29	13	27	10	24	08	22	05	19	02	16	30	14	28	12	25	09	23
AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT	AAI	ALDO-RIACT
BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT	ALDO-RIACT	CORT-CT2	BI-HNS-IRMA	CGA-RIACT
CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2	BI-HNS-IRMA	ELSA-ACTH	CEA-RIACT	CORT-CT2
CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH	CEA-RIACT	ELSA-CA125II	CGA-RIACT	ELSA-ACTH
CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II	CGA-RIACT	ELSA-PTH	CORT-CT2	ELSA-CA125II
ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH	CORT-CT2	ESTR-CTRIA	ELSA2-AFP	ELSA-PTH
ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA	ELSA2-AFP	ESTR-US-CT	ELSA-ACTH	ESTR-CTRIA
ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT	ELSA-ACTH	GASK-PR	ELSA-CA125II	ESTR-US-CT
ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR	ELSA-CA125II	HGH-RIACT	ELSA-CA15-3	GASK-PR
ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT	ELSA-CA15-3	IGF1-RIACT	ELSA-CA19-9	HGH-RIACT
ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT	ELSA-CA19-9	IRMA-HCT	ELSA-CA72-4	IGF1-RIACT
ELSA-CYFRA-	IRMA-CPEP	ELSA-CA72-4	OCFD03-FT4	ELSA-CYFRA-	IRMA-CPEP	ELSA-CA72-4	IRMA-HCT	ELSA-CYFRA-	IRMA-HCT	ELSA-CA72-4	IRMA-HCT	ELSA-CYFRA-	IRMA-CPEP	ELSA-CA72-4	OCFD03-FT4	ELSA-CYFRA-	IRMA-CPEP	ELSA-CA72-4	IRMA-HCT	ELSA-CYFRA-	IRMA-HCT	ELSA-CA72-4	IRMA-HCT	ELSA-CYFRA-	IRMA-CPEP
ELSA-FBHC	IRMA-HCT	ELSA-CYFRA-	OCFD03-FT4-K	ELSA-FBHC	IRMA-HCT	ELSA-CYFRA-	OCFD03-FT4	ELSA-FBHC	OCFA07-AFP	ELSA-CYFRA-	OCFD03-FT4	ELSA-FBHC	IRMA-HCT	ELSA-CYFRA-	OCFD03-FT4-K	ELSA-FBHC	IRMA-HCT	ELSA-CYFRA-	OCFD03-FT4	ELSA-FBHC	OCFA07-AFP	ELSA-CYFRA-	OCFD03-FT4	ELSA-FBHC	IRMA-HCT
ELSA-NSE	OCFA07-AFP	ELSA-FBHC	OCFD53-FT4	ELSA-NSE	OCFA07-AFP	ELSA-FBHC	OCFD03-FT4-K	ELSA-NSE	OCFD03-FT4	ELSA-NSE	OCFD03-FT4	ELSA-NSE	OCFA07-AFP	ELSA-FBHC	OCFD03-FT4	ELSA-NSE	OCFA07-AFP	ELSA-FBHC	OCFD03-FT4-K	ELSA-NSE	OCFD03-FT4	ELSA-FBHC	OCFD03-FT4-K	ELSA-NSE	OCFA07-AFP
ESTR-CTRIA	OCFD03-FT4	ELSA-NSE	OCFE07-TRYPS	ESTR-CTRIA	OCFD03-FT4	ELSA-NSE	OCFD03-FT4-K	ESTR-CTRIA	OCFD03-FT4	ELSA-NSE	OCFD03-FT4	ESTR-CTRIA	OCFD03-FT4	ELSA-NSE	OCFD03-FT4	ESTR-CTRIA	OCFD03-FT4	ELSA-NSE	OCFD03-FT4	ESTR-CTRIA	OCFD03-FT4-K	ELSA-NSE	OCFD03-FT4	ESTR-CTRIA	OCFD03-FT4
FPSA-RIACT	OCFD03-FT4-K	FPSA-RIACT	OCFH07-FT3	FPSA-RIACT	OCFD03-FT4-K	FPSA-RIACT	OCFE07-TRYPS	FPSA-RIACT	OCFD53-FT4	FPSA-RIACT	OCFD07-TRYPS	FPSA-RIACT	OCFD03-FT4-K	FPSA-RIACT	OCFH07-FT3	FPSA-RIACT	OCFD03-FT4-K	FPSA-RIACT	OCFE07-TRYPS	FPSA-RIACT	OCFD53-FT4	FPSA-RIACT	OCFE07-TRYPS	FPSA-RIACT	OCFD03-FT4-K
HGH-RIACT	OCFD53-FT4	HGH-RIACT	OCFH57-FT3	HGH-RIACT	OCFD53-FT4	HGH-RIACT	OCFH07-FT3	HGH-RIACT	OCFE07-TRYPS	HGH-RIACT	OCFH07-FT3	HGH-RIACT	OCFD53-FT4	HGH-RIACT	OCFH57-FT3	HGH-RIACT	OCFD53-FT4	HGH-RIACT	OCFH07-FT3	HGH-RIACT	OCFE07-TRYPS	HGH-RIACT	OCFH07-FT3	HGH-RIACT	OCFD53-FT4
IRMA-HCT	OCFE07-TRYPS	IGF1-RIACT	OCPE07-T3	IRMA-CPEP	OCFE07-TRYPS	IGF1-RIACT	OCFH57-FT3	IRMA-CPEP	OCFH07-FT3	IGF1-RIACT	OCFH57-FT3	IRMA-HCT	OCFE07-TRYPS	IGF1-RIACT	OCPE07-T3	IRMA-CPEP	OCFE07-TRYPS	IGF1-RIACT	OCFH57-FT3	IRMA-CPEP	OCFH07-FT3	IGF1-RIACT	OCFH57-FT3	IRMA-HCT	OCFE07-TRYPS
OCFA07-AFP	OCFH07-FT3	IRMA-CPEP	OCPE57-T3	IRMA-HCT	OCFH07-FT3	IRMA-HCT	OCFM07-CA50	IRMA-HCT	OCFH57-FT3	IRMA-CPEP	OCFM07-CA50	OCFA07-AFP	OCFH07-FT3	IRMA-CPEP	OCPE57-T3	IRMA-HCT	OCFH07-FT3	IRMA-HCT	OCFM07-CA50	IRMA-HCT	OCFH57-FT3	IRMA-CPEP	OCFM07-CA50	OCFA07-AFP	OCFH07-FT3
OCFD03-FT4	OCFH57-FT3	IRMA-HCT	OCPG07-T4	OCFA07-AFP	OCFH57-FT3	OCFA07-AFP	OCPE07-T3	OCFA07-AFP	OCPE07-T3	IRMA-HCT	OCPE07-T3	OCFD03-FT4	OCFH57-FT3	IRMA-HCT	OCPG07-T4	OCFA07-AFP	OCFH57-FT3	OCFA07-AFP	OCPE07-T3	OCFA07-AFP	OCPE07-T3	OCFA07-AFP	OCPE07-T3	OCFA07-AFP	OCPE07-T3
OCFD03-FT4-K	OCFM07-CA50	OCFA07-AFP	OCPG57-T4	OCFD03-FT4	OCFM07-CA50	OCFD03-FT4	OCPE57-T3	OCFD03-FT4	OCPE57-T3	OCFA07-AFP	OCPE57-T3	OCFD03-FT4-K	OCFM07-CA50	OCFA07-AFP	OCPG57-T4	OCFD03-FT4	OCFM07-CA50	OCFD03-FT4	OCPE57-T3	OCFD03-FT4	OCPE57-T3	OCFA07-AFP	OCPE57-T3	OCFD03-FT4-K	OCFM07-CA50
OCFD53-FT4	OCPE07-T3	OCFD03-FT4	OCPL07-TSH	OCFD03-FT4-K	OCPE07-T3	OCFD03-FT4-K	OCPG07-T4	OCFD03-FT4-K	OCPG07-T4	OCFD03-FT4	OCPE07-T3	OCFD03-FT4	OCPL07-TSH	OCFD03-FT4-K	OCPE07-T3	OCFD03-FT4	OCPL07-TSH	OCFD03-FT4-K	OCPE07-T3	OCFD03-FT4	OCPE07-T3	OCFA07-AFP	OCPE57-T3	OCFD03-FT4-K	OCFM07-CA50
OCFE07-TRYPS	OCPE57-T3	OCFD03-FT4-K	OCPL57-TSH	OCFD53-FT4	OCPE57-T3	OCFD53-FT4	OCPG57-T4	OCFD53-FT4	OCPE57-T4	OCFD53-FT4	OCPE57-T3	OCFD03-FT4-K	OCPL57-TSH	OCFD53-FT4	OCPE57-T3	OCFD53-FT4	OCPL57-TSH	OCFD53-FT4	OCPE57-T3	OCFD53-FT4	OCPE57-T4	OCFD03-FT4-K	OCPE57-T4	OCFD03-FT4-K	OCPE57-T3
OCFH07-FT3	OCPG07-T4	OCFD53-FT4	OCPM03-HCG	OCFE07-TRYPS	OCPG07-T4	OCFE07-TRYPS	OCPL07-TSH	OCFE07-TRYPS	OCPL07-TSH	OCFD53-FT4	OCPL07-TSH	OCFH07-FT3	OCPG07-T4	OCFD53-FT4	OCPM03-HCG	OCFE07-TRYPS	OCPG07-T4	OCFE07-TRYPS	OCPL07-TSH	OCFE07-TRYPS	OCPL07-TSH	OCFD53-FT4	OCPL07-TSH	OCFH07-FT3	OCPG07-T4
OCFH57-FT3	OCPG57-T4	OCFE07-TRYPS	OCRU03-TBG	OCFH07-FT3	OCPG57-T4	OCFH07-FT3	OCPL57-TSH	OCFH07-FT3	OCPL57-TSH	OCFE07-TRYPS	OCPL57-TSH	OCFH57-FT3	OCPG57-T4	OCFE07-TRYPS	OCRU03-TBG	OCFH07-FT3	OCPG57-T4	OCFH07-FT3	OCPL57-TSH	OCFH07-FT3	OCPL57-TSH	OCFE07-TRYPS	OCPL57-TSH	OCFH57-FT3	OCPG57-T4
OCFK07-PIIP	OCPL07-TSH	OCFH07-FT3	OSTEO-RIACT	OCFH57-FT3	OCPL07-TSH	OCFH57-FT3	OCPM03-HCG	OCFH57-FT3	OCPM03-HCG	OCFH07-FT3	OCPM03-HCG	OCFK07-PIIP	OCPL07-TSH	OCFH07-FT3	OSTEO-RIACT	OCFH57-FT3	OCPL07-TSH	OCFH57-FT3	OCPM03-HCG	OCFH57-FT3	OCPM03-HCG	OCFH07-FT3	OCPL07-TSH	OCFH57-FT3	OCPG57-T4
OCPE07-T3	OCPL57-TSH	OCFH57-FT3	RENINE	OCFK07-PIIP	OCPL57-TSH	OCFK07-PIIP	OCRU03-TBG	OCFK07-PIIP	OCRU03-TBG	OCFH57-FT3	OCRU03-TBG	OCPE07-T3	OCPL57-TSH	OCFH57-FT3	RENINE	OCFK07-PIIP	OCPL57-TSH	OCFK07-PIIP	OCRU03-TBG	OCFK07-PIIP	OCRU03-TBG	OCFH57-FT3	OCRU03-TBG	OCPE07-T3	OCPL57-TSH
OCPE57-T3	OCPM03-HCG	OCFK07-PIIP	R-GM-100	OCFM07-CA50	OCPM03-HCG	OCPE07-T3	OSTEO-RIACT	OCFM07-CA50	OSTEO-RIACT	OCFK07-PIIP	OSTEO-RIACT	OCPE57-T3	OCPM03-HCG	OCFK07-PIIP	R-GM-100	OCFM07-CA50	OCPM03-HCG	OCPE07-T3	OSTEO-RIACT	OCFM07-CA50	OSTEO-RIACT	OCFK07-PIIP	OSTEO-RIACT	OCPE57-T3	OCPM03-HCG
OCPG07-T4	OCRU03-TBG	OCFM07-CA50	SHBG-RIACT	OCPE07-T3	OCRU03-TBG	OCPE57-T3	RENINE	OCPE07-T3	RENINE	OCFM07-CA50	RENINE	OCPG07-T4	OCRU03-TBG	OCFM07-CA50	SHBG-RIACT	OCPE07-T3	OCRU03-TBG	OCPE57-T3	RENINE	OCPE07-T3	RENINE	OCFM07-CA50	RENINE	OCPG07-T4	OCRU03-TBG
OCPG57-T4	OSTEO-RIACT	OCPE07-T3	TESTO-CT2	OCPE57-T3	OSTEO-RIACT	OCPG07-T4	R-GM-100	OCPE57-T3	R-GM-100	OCPE07-T3	R-GM-100	OCPG57-T4	OSTEO-RIACT	OCPE07-T3	TESTO-CT2	OCPE57-T3	OSTEO-RIACT	OCPG07-T4	R-GM-100	OCPE57-T3	TESTO-CT2	OCPE57-T3	R-GM-100	OCPG57-T4	OSTEO-RIACT
OCPL07-TSH	RENINE	OCPE57-T3	THYRO	OCPE07-T3	THYRO	OCPE57-T4	SHBG-RIACT	OCPE07-T4	SHBG-RIACT	OCPE07-T3	SHBG-RIACT	OCPL07-TSH	RENINE	OCPE57-T3	THYRO	OCPE07-T4	SHBG-RIACT	OCPE57-T3	THYRO	OCPE07-T4	SHBG-RIACT	OCPE57-T3	THYRO	OCPE07-T4	SHBG-RIACT
OCPL57-TSH	R-GM-100	OCPE07-T4		OCPE57-T4	R-GM-100	OCPL07-TSH	TESTO-CT2	OCPE57-T4	TESTO-CT2	OCPE07-T4	TESTO-CT2	OCPL57-TSH	R-GM-100	OCPE07-T4		OCPE57-T4	R-GM-100	OCPL07-TSH	TESTO-CT2	OCPE57-T4	TESTO-CT2	OCPE07-T4	TESTO-CT2	OCPL57-TSH	R-GM-100
OCPM03-HCG	SHBG-RIACT	OCPE57-T4		OCPL07-TSH	SHBG-RIACT	OCPL57-TSH	THYRO	OCPL07-TSH	THYRO	OCPE57-T4	THYRO	OCPM03-HCG	SHBG-RIACT	OCPE57-T4		OCPL07-TSH	SHBG-RIACT	OCPL57-TSH	THYRO	OCPL07-TSH	THYRO	OCPE57-T4	THYRO	OCPM03-HCG	SHBG-RIACT
OCRU03-TBG	TESTO-CT2	OCPL07-TSH		OCPL57-TSH	TESTO-CT2	OCRU03-TBG		OCPL57-TSH		OCPL07-TSH		OCRU03-TBG	TESTO-CT2	OCPL07-TSH		OCPL57-TSH	TESTO-CT2	OCRU03-TBG		OCPL57-TSH		OCPL07-TSH		OCRU03-TBG	TESTO-CT2
PSA-RIACT	THYRO	OCPL57-TSH		OCPM03-HCG	THYRO	PSA-RIACT		OCPM03-HCG		OCPL57-TSH		PSA-RIACT	THYRO	OCPL57-TSH		OCPM03-HCG	THYRO	PSA-RIACT		OCPM03-HCG		OCPL57-TSH		PSA-RIACT	THYRO
R-GM-100		OCRU03-TBG		OCRU03-TBG		R-GM-100		OCRU03-TBG		R-GM-100		R-GM-100		OCRU03-TBG		R-GM-100		OCRU03-TBG		R-GM-100		OCRU03-TBG		R-GM-100	
SHBG-RIACT		PSA-RIACT		PSA-RIACT		TESTO-CT2		PSA-RIACT		SHBG-RIACT		SHBG-RIACT		PSA-RIACT		SHBG-RIACT		PSA-RIACT		SHBG-RIACT		PSA-RIACT		SHBG-RIACT	
TESTO-CT2		R-GM-100		R-GM-100		TESTO-CT2		R-GM-100		TESTO-CT2		TESTO-CT2		R-GM-100		TESTO-CT2		R-GM-100		TESTO-CT2		R-GM-100		TESTO-CT2	
		TESTO-CT2		SHBG-RIACT		SHBG-RIACT		SHBG-RIACT		TESTO-CT2		SHBG-RIACT		SHBG-RIACT		TESTO-CT2		SHBG-RIACT		SHBG-RIACT		TESTO-CT2		SHBG-RIACT	
		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2		TESTO-CT2	



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