

CALENDRIER DE COMMANDE

2022

PRODUITS RIA

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
43	CGA-RIACT	22/10/2021	10/11/2021
44	OCFA07-AFP	26/10/2021	10/11/2021
44	OCRU03-TBG	29/10/2021	17/11/2021
45	ALDO-RIACT	05/11/2021	22/11/2021
45	ALDO-US	05/11/2021	22/11/2021
44	IGF1-RIACT	29/10/2021	22/11/2021
44	IGF1-US	29/10/2021	22/11/2021
46	CA50-US	08/11/2021	24/11/2021
45	CPEP-US	05/11/2021	24/11/2021
45	IRMA-CPEP	05/11/2021	24/11/2021
46	OCFM07-CA50	08/11/2021	24/11/2021
46	ELSA2-AFP	12/11/2021	01/12/2021
46	ELSA-CA15-3	10/11/2021	01/12/2021
46	ELSA-CA19-9	10/11/2021	01/12/2021
46	ELSA-CA72-4	10/11/2021	01/12/2021
46	ELSA-FBHCG	10/11/2021	01/12/2021
45	MSK/TR	04/11/2021	01/12/2021
47	OCFE07-TRYPS	15/11/2021	01/12/2021
45	RBAN/25	04/11/2021	01/12/2021
44	TR-AB-CT	27/10/2021	01/12/2021
47	TRYPS-DIA	15/11/2021	01/12/2021
47	TRYPS-US	15/11/2021	01/12/2021
47	BI-INS-IRMA	15/11/2021	03/12/2021
47	BI-INSU-US	15/11/2021	03/12/2021
46	ELSA-CYFRA-	10/11/2021	03/12/2021
47	ELSA-NSE	16/11/2021	03/12/2021
47	OCFK07-PIIIP	17/11/2021	03/12/2021
47	PIIIP-US	17/11/2021	03/12/2021
47	AAI	16/11/2021	06/12/2021
47	AAI-US	16/11/2021	06/12/2021
47	CEA-RIACT	15/11/2021	06/12/2021
46	ACHR-AB	10/11/2021	08/12/2021
47	ELSA-CA125II	17/11/2021	08/12/2021
48	FPSA-RIACT	22/11/2021	08/12/2021
48	IRMA-HCT	22/11/2021	08/12/2021
48	PSA-RIACT	22/11/2021	08/12/2021
46	TPO-AB-CT	10/11/2021	08/12/2021
48	OCFD03-FT4	24/11/2021	10/12/2021
48	OCFD03-FT4-K	24/11/2021	10/12/2021
48	OCFD53-FT4	24/11/2021	10/12/2021
48	OCFH07-FT3	24/11/2021	10/12/2021
48	OCFH57-FT3	24/11/2021	10/12/2021
48	OCPL07-TSH	23/11/2021	10/12/2021
48	OCPL57-TSH	23/11/2021	10/12/2021
48	CORT-CT2	25/11/2021	15/12/2021
48	CORT-US	25/11/2021	15/12/2021

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
49	GASK-PR	29/11/2021	15/12/2021
49	GASK-PR-US	29/11/2021	15/12/2021
49	OCPE07-T3	30/11/2021	15/12/2021
49	OCPE57-T3	30/11/2021	15/12/2021
49	OCPG07-T4	30/11/2021	15/12/2021
49	OCPG57-T4	30/11/2021	15/12/2021
49	R-GM-100	29/11/2021	15/12/2021
49	T3-US	30/11/2021	15/12/2021
49	T4-US	30/11/2021	15/12/2021
47	TGAB I STEP	15/11/2021	16/12/2021
48	ESTR-S-US	25/11/2021	17/12/2021
48	ESTR-US-CT	25/11/2021	17/12/2021
49	OSTEO-RIACT	02/12/2021	17/12/2021
49	OSTEO-US	02/12/2021	17/12/2021
49	SHBG-RIACT	29/11/2021	17/12/2021
48	TESTO-CT2	25/11/2021	17/12/2021
48	TESTO-US	25/11/2021	17/12/2021
49	ESTR-CTRIA	02/12/2021	20/12/2021
49	ESTR-US	02/12/2021	20/12/2021
49	RENINE	30/11/2021	20/12/2021
49	ELSA-ACTH	01/12/2021	22/12/2021
50	ELSA-PTH	06/12/2021	22/12/2021
49	HCG-US	03/12/2021	22/12/2021
49	HGH-RIACT	03/12/2021	22/12/2021
49	HGH-US	03/12/2021	22/12/2021
50	IRMA-HCT	06/12/2021	22/12/2021
48	MSK/TR	25/11/2021	22/12/2021
49	OCPM03-HCG	03/12/2021	22/12/2021
48	RBAN/25	25/11/2021	22/12/2021
50	THYRO	06/12/2021	22/12/2021
47	TR-AB-CT	18/11/2021	22/12/2021
50	ELSA-FBHCGSP	09/12/2021	24/12/2021
50	ELSA2-AFP	10/12/2021	29/12/2021
50	ELSA-CA15-3	08/12/2021	29/12/2021
50	ELSA-CA19-9	08/12/2021	29/12/2021
50	ELSA-CA72-4	09/12/2021	29/12/2021
50	ELSA-FBHCG	09/12/2021	29/12/2021
51	OCFE07-TRYPS	13/12/2021	29/12/2021
51	TRYPS-DIA	13/12/2021	29/12/2021
51	TRYPS-US	13/12/2021	29/12/2021
49	17-OHP-CT	29/11/2021	30/12/2021
49	PROGEST-RIA	29/11/2021	30/12/2021
49	REN-CT2P	29/11/2021	30/12/2021
51	BI-INS-IRMA	13/12/2021	31/12/2021
51	BI-INSU-US	13/12/2021	31/12/2021
50	ELSA-CYFRA-	09/12/2021	31/12/2021

CALENDRIER DE COMMANDE

2022

PRODUITS RIA

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
51	ELSA-NSE	14/12/2021	31/12/2021
51	OCFK07-PIIIP	15/12/2021	31/12/2021
51	PIIIP-US	15/12/2021	31/12/2021
51	AAI	14/12/2021	03/01/2022
51	AAI-US	14/12/2021	03/01/2022
51	CEA-RIACT	13/12/2021	03/01/2022
52	CA50-US	20/12/2021	05/01/2022
51	CGA-RIACT	17/12/2021	05/01/2022
51	CPEP-US	16/12/2021	05/01/2022
51	ELSA-CA125II	15/12/2021	05/01/2022
52	FPSA-RIACT	20/12/2021	05/01/2022
51	IRMA-CPEP	16/12/2021	05/01/2022
52	IRMA-HCT	20/12/2021	05/01/2022
52	OCFA07-AFP	21/12/2021	05/01/2022
52	OCFM07-CA50	20/12/2021	05/01/2022
52	PSA-RIACT	20/12/2021	05/01/2022
52	OCFD03-FT4	22/12/2021	07/01/2022
52	OCFD03-FT4-K	22/12/2021	07/01/2022
52	OCFD53-FT4	22/12/2021	07/01/2022
52	OCFH07-FT3	22/12/2021	07/01/2022
52	OCFH57-FT3	22/12/2021	07/01/2022
52	OCPL07-TSH	21/12/2021	07/01/2022
52	OCPL57-TSH	21/12/2021	07/01/2022
49	ACHR-AB	01/12/2021	12/01/2022
52	CORT-CT2	23/12/2021	12/01/2022
52	CORT-US	23/12/2021	12/01/2022
53	GASK-PR	27/12/2021	12/01/2022
53	GASK-PR-US	27/12/2021	12/01/2022
53	OCPE07-T3	28/12/2021	12/01/2022
53	OCPE57-T3	28/12/2021	12/01/2022
53	OCPG07-T4	28/12/2021	12/01/2022
53	OCPG57-T4	28/12/2021	12/01/2022
53	OCRU03-TBG	27/12/2021	12/01/2022
53	R-GM-100	27/12/2021	12/01/2022
53	T3-US	28/12/2021	12/01/2022
53	T4-US	28/12/2021	12/01/2022
51	TGAB I STEP	13/12/2021	13/01/2022
52	ESTR-S-US	23/12/2021	14/01/2022
52	ESTR-US-CT	23/12/2021	14/01/2022
53	OSTEO-RIACT	30/12/2021	14/01/2022
53	OSTEO-US	30/12/2021	14/01/2022
53	ALDO-RIACT	30/12/2021	17/01/2022
53	ALDO-US	30/12/2021	17/01/2022
53	IGF1-RIACT	27/12/2021	17/01/2022
53	IGF1-US	27/12/2021	17/01/2022
53	RENINE	28/12/2021	17/01/2022
53	ELSA-ACTH	29/12/2021	19/01/2022
2	ELSA-PTH	03/01/2022	19/01/2022
53	HGH-RIACT	31/12/2021	19/01/2022
53	HGH-US	31/12/2021	19/01/2022
2	IRMA-HCT	03/01/2022	19/01/2022
2	THYRO	03/01/2022	19/01/2022
2	ELSA-FBHCGSP	06/01/2022	21/01/2022
2	ELSA2-AFP	07/01/2022	26/01/2022
2	ELSA-CA15-3	05/01/2022	26/01/2022

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
2	ELSA-CA19-9	05/01/2022	26/01/2022
2	ELSA-CA72-4	06/01/2022	26/01/2022
2	ELSA-FBHCG	06/01/2022	26/01/2022
53	MSK/TR	30/12/2021	26/01/2022
3	OCFE07-TRYPS	10/01/2022	26/01/2022
53	RBAN/25	30/12/2021	26/01/2022
52	TR-AB-CT	23/12/2021	26/01/2022
3	TRYPS-DIA	10/01/2022	26/01/2022
3	TRYPS-US	10/01/2022	26/01/2022
51	VGC/25	15/12/2021	26/01/2022
51	VGK/25	15/12/2021	26/01/2022
53	REN-CT2P	27/12/2021	27/01/2022
3	BI-INS-IRMA	10/01/2022	28/01/2022
3	BI-INSU-US	10/01/2022	28/01/2022
2	ELSA-CYFRA-	06/01/2022	28/01/2022
3	ELSA-NSE	11/01/2022	28/01/2022
3	OCFK07-PIIIP	12/01/2022	28/01/2022
3	PIIIP-US	12/01/2022	28/01/2022
3	AAI	11/01/2022	31/01/2022
3	AAI-US	11/01/2022	31/01/2022
3	CEA-RIACT	10/01/2022	31/01/2022
2	ACHR-AB	05/01/2022	02/02/2022
3	ELSA-CA125II	12/01/2022	02/02/2022
4	FPSA-RIACT	17/01/2022	02/02/2022
4	IRMA-HCT	17/01/2022	02/02/2022
4	PSA-RIACT	17/01/2022	02/02/2022
2	TPO-AB-CT	05/01/2022	02/02/2022
4	OCFD03-FT4	19/01/2022	04/02/2022
4	OCFD03-FT4-K	19/01/2022	04/02/2022
4	OCFD53-FT4	19/01/2022	04/02/2022
4	OCFH07-FT3	19/01/2022	04/02/2022
4	OCFH57-FT3	19/01/2022	04/02/2022
4	OCPL07-TSH	18/01/2022	04/02/2022
4	OCPL57-TSH	18/01/2022	04/02/2022
4	CORT-CT2	20/01/2022	09/02/2022
4	CORT-US	20/01/2022	09/02/2022
5	GASK-PR	24/01/2022	09/02/2022
5	GASK-PR-US	24/01/2022	09/02/2022
5	OCPE07-T3	25/01/2022	09/02/2022
5	OCPE57-T3	25/01/2022	09/02/2022
5	OCPG07-T4	25/01/2022	09/02/2022
5	OCPG57-T4	25/01/2022	09/02/2022
5	R-GM-100	24/01/2022	09/02/2022
5	T3-US	25/01/2022	09/02/2022
5	T4-US	25/01/2022	09/02/2022
3	17-OHP-CT	10/01/2022	10/02/2022
3	PROGEST-RIA	10/01/2022	10/02/2022
3	TGAB I STEP	10/01/2022	10/02/2022
4	ESTR-S-US	20/01/2022	11/02/2022
4	ESTR-US-CT	20/01/2022	11/02/2022

CALENDRIER DE COMMANDE

2022

PRODUITS RIA

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
5	OSTEO-RIACT	27/01/2022	11/02/2022
5	OSTEO-US	27/01/2022	11/02/2022
5	SHBG-RIACT	24/01/2022	11/02/2022
4	TESTO-CT2	20/01/2022	11/02/2022
4	TESTO-US	20/01/2022	11/02/2022
5	ESTR-CTRIA	27/01/2022	14/02/2022
5	ESTR-US	27/01/2022	14/02/2022
5	RENINE	25/01/2022	14/02/2022
6	CA50-US	31/01/2022	16/02/2022
5	CPEP-US	27/01/2022	16/02/2022
5	ELSA-ACTH	26/01/2022	16/02/2022
6	ELSA-PTH	31/01/2022	16/02/2022
5	HCG-US	28/01/2022	16/02/2022
5	HGH-RIACT	28/01/2022	16/02/2022
5	HGH-US	28/01/2022	16/02/2022
5	IRMA-CPEP	27/01/2022	16/02/2022
6	IRMA-HCT	31/01/2022	16/02/2022
6	OCFM07-CA50	31/01/2022	16/02/2022
5	OCPM03-HCG	28/01/2022	16/02/2022
6	THYRO	31/01/2022	16/02/2022
6	ELSA-FBHCGSP	03/02/2022	18/02/2022
6	ELSA2-AFP	04/02/2022	23/02/2022
6	ELSA-CA15-3	02/02/2022	23/02/2022
6	ELSA-CA19-9	02/02/2022	23/02/2022
6	ELSA-CA72-4	03/02/2022	23/02/2022
6	ELSA-FBHCG	03/02/2022	23/02/2022
5	MSK/TR	27/01/2022	23/02/2022
7	OCFE07-TRYPS	07/02/2022	23/02/2022
5	RBAN/25	27/01/2022	23/02/2022
4	TR-AB-CT	20/01/2022	23/02/2022
7	TRYPS-DIA	07/02/2022	23/02/2022
7	TRYPS-US	07/02/2022	23/02/2022
5	REN-CT2P	24/01/2022	24/02/2022
7	BI-INS-IRMA	07/02/2022	25/02/2022
7	BI-INSU-US	07/02/2022	25/02/2022
6	ELSA-CYFRA-	03/02/2022	25/02/2022
7	ELSA-NSE	08/02/2022	25/02/2022
7	OCFK07-PIIIP	09/02/2022	25/02/2022
7	PIIIP-US	09/02/2022	25/02/2022
7	AAI	08/02/2022	28/02/2022
7	AAI-US	08/02/2022	28/02/2022
7	CEA-RIACT	07/02/2022	28/02/2022
6	ACHR-AB	02/02/2022	02/03/2022
7	CGA-RIACT	11/02/2022	02/03/2022
7	ELSA-CA125II	09/02/2022	02/03/2022
8	FPSA-RIACT	14/02/2022	02/03/2022
8	IRMA-HCT	14/02/2022	02/03/2022
8	OCFA07-AFP	15/02/2022	02/03/2022
8	PSA-RIACT	14/02/2022	02/03/2022
8	OCFD03-FT4	16/02/2022	04/03/2022
8	OCFD03-FT4-K	16/02/2022	04/03/2022
8	OCFD53-FT4	16/02/2022	04/03/2022
8	OCFH07-FT3	16/02/2022	04/03/2022
8	OCFH57-FT3	16/02/2022	04/03/2022
8	OCPL07-TSH	15/02/2022	04/03/2022

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
8	OCPL57-TSH	15/02/2022	04/03/2022
8	CORT-CT2	17/02/2022	09/03/2022
8	CORT-US	17/02/2022	09/03/2022
9	GASK-PR	21/02/2022	09/03/2022
9	GASK-PR-US	21/02/2022	09/03/2022
9	OCPE07-T3	22/02/2022	09/03/2022
9	OCPE57-T3	22/02/2022	09/03/2022
9	OCPG07-T4	22/02/2022	09/03/2022
9	OCPG57-T4	22/02/2022	09/03/2022
9	OCRU03-TBG	21/02/2022	09/03/2022
9	R-GM-100	21/02/2022	09/03/2022
9	T3-US	22/02/2022	09/03/2022
9	T4-US	22/02/2022	09/03/2022
7	VGC/25	09/02/2022	09/03/2022
7	VGK/25	09/02/2022	09/03/2022
7	TGAB I STEP	07/02/2022	10/03/2022
8	ESTR-S-US	17/02/2022	11/03/2022
8	ESTR-US-CT	17/02/2022	11/03/2022
9	OSTEO-RIACT	24/02/2022	11/03/2022
9	OSTEO-US	24/02/2022	11/03/2022
9	ALDO-RIACT	24/02/2022	14/03/2022
9	ALDO-US	24/02/2022	14/03/2022
9	IGF1-RIACT	21/02/2022	14/03/2022
9	IGF1-US	21/02/2022	14/03/2022
9	RENINE	22/02/2022	14/03/2022
9	ELSA-ACTH	23/02/2022	16/03/2022
10	ELSA-PTH	28/02/2022	16/03/2022
9	HGH-RIACT	25/02/2022	16/03/2022
9	HGH-US	25/02/2022	16/03/2022
10	IRMA-HCT	28/02/2022	16/03/2022
10	THYRO	28/02/2022	16/03/2022
10	ELSA-FBHCGSP	03/03/2022	18/03/2022
10	ELSA2-AFP	04/03/2022	23/03/2022
10	ELSA-CA15-3	02/03/2022	23/03/2022
10	ELSA-CA19-9	02/03/2022	23/03/2022
10	ELSA-CA72-4	03/03/2022	23/03/2022
10	ELSA-FBHCG	03/03/2022	23/03/2022
9	MSK/TR	24/02/2022	23/03/2022
11	OCFE07-TRYPS	07/03/2022	23/03/2022
9	RBAN/25	24/02/2022	23/03/2022
8	TR-AB-CT	17/02/2022	23/03/2022
11	TRYPS-DIA	07/03/2022	23/03/2022
11	TRYPS-US	07/03/2022	23/03/2022
9	17-OHP-CT	21/02/2022	24/03/2022
9	PROGEST-RIA	21/02/2022	24/03/2022
9	REN-CT2P	21/02/2022	24/03/2022
11	BI-INS-IRMA	07/03/2022	25/03/2022
11	BI-INSU-US	07/03/2022	25/03/2022
10	ELSA-CYFRA-	03/03/2022	25/03/2022
11	ELSA-NSE	08/03/2022	25/03/2022

CALENDRIER DE COMMANDE

2022

PRODUITS RIA

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
11	OCFK07-PIIIP	09/03/2022	25/03/2022
11	PIIIP-US	09/03/2022	25/03/2022
11	AAI	08/03/2022	28/03/2022
11	AAI-US	08/03/2022	28/03/2022
11	CEA-RIACT	07/03/2022	28/03/2022
10	ACHR-AB	02/03/2022	30/03/2022
12	CA50-US	14/03/2022	30/03/2022
11	CPEP-US	10/03/2022	30/03/2022
11	ELSA-CA125II	09/03/2022	30/03/2022
12	FPSA-RIACT	14/03/2022	30/03/2022
11	IRMA-CPEP	10/03/2022	30/03/2022
12	IRMA-HCT	14/03/2022	30/03/2022
12	OCFM07-CA50	14/03/2022	30/03/2022
12	PSA-RIACT	14/03/2022	30/03/2022
10	TPO-AB-CT	02/03/2022	30/03/2022
12	OCFD03-FT4	16/03/2022	01/04/2022
12	OCFD03-FT4-K	16/03/2022	01/04/2022
12	OCFD53-FT4	16/03/2022	01/04/2022
12	OCFH07-FT3	16/03/2022	01/04/2022
12	OCFH57-FT3	16/03/2022	01/04/2022
12	OCPL07-TSH	15/03/2022	01/04/2022
12	OCPL57-TSH	15/03/2022	01/04/2022
12	CORT-CT2	17/03/2022	06/04/2022
12	CORT-US	17/03/2022	06/04/2022
13	GASK-PR	21/03/2022	06/04/2022
13	GASK-PR-US	21/03/2022	06/04/2022
13	OCPE07-T3	22/03/2022	06/04/2022
13	OCPE57-T3	22/03/2022	06/04/2022
13	OCPG07-T4	22/03/2022	06/04/2022
13	OCPG57-T4	22/03/2022	06/04/2022
13	R-GM-100	21/03/2022	06/04/2022
13	T3-US	22/03/2022	06/04/2022
13	T4-US	22/03/2022	06/04/2022
11	TGAB I STEP	07/03/2022	07/04/2022
12	ESTR-S-US	17/03/2022	08/04/2022
12	ESTR-US-CT	17/03/2022	08/04/2022
13	OSTEO-RIACT	24/03/2022	08/04/2022
13	OSTEO-US	24/03/2022	08/04/2022
13	SHBG-RIACT	21/03/2022	08/04/2022
12	TESTO-CT2	17/03/2022	08/04/2022
12	TESTO-US	17/03/2022	08/04/2022
13	ESTR-CTRIA	24/03/2022	11/04/2022
13	ESTR-US	24/03/2022	11/04/2022
13	RENINE	22/03/2022	11/04/2022
13	ELSA-ACTH	23/03/2022	13/04/2022
14	ELSA-PTH	28/03/2022	13/04/2022
13	HCG-US	25/03/2022	13/04/2022
13	HGH-RIACT	25/03/2022	13/04/2022
13	HGH-US	25/03/2022	13/04/2022
14	IRMA-HCT	28/03/2022	13/04/2022
13	OCPM03-HCG	25/03/2022	13/04/2022
14	THYRO	28/03/2022	13/04/2022
14	ELSA-FBHCGSP	31/03/2022	15/04/2022
13	MSK/TR	24/03/2022	20/04/2022
13	RBAN/25	24/03/2022	20/04/2022

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
12	TR-AB-CT	17/03/2022	20/04/2022
14	ELSA2-AFP	01/04/2022	21/04/2022
14	ELSA-CA15-3	30/03/2022	21/04/2022
14	ELSA-CA19-9	30/03/2022	21/04/2022
14	ELSA-CA72-4	31/03/2022	21/04/2022
14	ELSA-FBHCG	31/03/2022	21/04/2022
15	OCFE07-TRYPS	04/04/2022	21/04/2022
15	TRYPS-DIA	04/04/2022	21/04/2022
15	TRYPS-US	04/04/2022	21/04/2022
15	BI-INS-IRMA	04/04/2022	22/04/2022
15	BI-INSU-US	04/04/2022	22/04/2022
14	ELSA-CYFRA-	31/03/2022	22/04/2022
15	ELSA-NSE	05/04/2022	22/04/2022
15	OCFK07-PIIIP	06/04/2022	22/04/2022
15	PIIIP-US	06/04/2022	22/04/2022
13	REN-CT2P	21/03/2022	22/04/2022
14	AAI	02/04/2022	25/04/2022
14	AAI-US	02/04/2022	25/04/2022
15	CEA-RIACT	04/04/2022	25/04/2022
14	ACHR-AB	30/03/2022	27/04/2022
15	CGA-RIACT	08/04/2022	27/04/2022
15	ELSA-CA125II	06/04/2022	27/04/2022
16	FPSA-RIACT	12/04/2022	27/04/2022
16	IRMA-HCT	12/04/2022	27/04/2022
16	OCFA07-AFP	12/04/2022	27/04/2022
16	PSA-RIACT	12/04/2022	27/04/2022
16	OCFD03-FT4	13/04/2022	29/04/2022
16	OCFD03-FT4-K	13/04/2022	29/04/2022
16	OCFD53-FT4	13/04/2022	29/04/2022
16	OCFH07-FT3	13/04/2022	29/04/2022
16	OCFH57-FT3	13/04/2022	29/04/2022
16	OCPL07-TSH	12/04/2022	29/04/2022
16	OCPL57-TSH	12/04/2022	29/04/2022
16	CORT-CT2	14/04/2022	04/05/2022
16	CORT-US	14/04/2022	04/05/2022
17	GASK-PR	18/04/2022	04/05/2022
17	GASK-PR-US	18/04/2022	04/05/2022
17	OCPE07-T3	19/04/2022	04/05/2022
17	OCPE57-T3	19/04/2022	04/05/2022
17	OCPG07-T4	19/04/2022	04/05/2022
17	OCPG57-T4	19/04/2022	04/05/2022
16	OCRU03-TBG	15/04/2022	04/05/2022
16	R-GM-100	15/04/2022	04/05/2022
17	T3-US	19/04/2022	04/05/2022
17	T4-US	19/04/2022	04/05/2022
15	VGC/25	06/04/2022	04/05/2022
15	VGK/25	06/04/2022	04/05/2022
15	17-OHP-CT	05/04/2022	05/05/2022
15	PROGEST-RIA	05/04/2022	05/05/2022
15	TGAB I STEP	04/04/2022	05/05/2022

CALENDRIER DE COMMANDE

2022


PRODUITS RIA

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
16	ESTR-S-US	14/04/2022	06/05/2022
16	ESTR-US-CT	14/04/2022	06/05/2022
17	ALDO-RIACT	21/04/2022	09/05/2022
17	ALDO-US	21/04/2022	09/05/2022
16	IGF1-RIACT	15/04/2022	09/05/2022
16	IGF1-US	15/04/2022	09/05/2022
17	RENINE	19/04/2022	09/05/2022
18	CA50-US	25/04/2022	11/05/2022
17	CPEP-US	21/04/2022	11/05/2022
17	ELSA-ACTH	20/04/2022	11/05/2022
18	ELSA-PTH	25/04/2022	11/05/2022
17	ESTR-CTRIA	21/04/2022	11/05/2022
17	ESTR-US	21/04/2022	11/05/2022
17	HCG-US	22/04/2022	11/05/2022
17	HGH-RIACT	22/04/2022	11/05/2022
17	HGH-US	22/04/2022	11/05/2022
17	IRMA-CPEP	21/04/2022	11/05/2022
18	IRMA-HCT	25/04/2022	11/05/2022
18	OCFM07-CA50	25/04/2022	11/05/2022
17	OCPM03-HCG	22/04/2022	11/05/2022
18	OSTEO-RIACT	28/04/2022	13/05/2022
18	OSTEO-US	28/04/2022	13/05/2022
18	SHBG-RIACT	25/04/2022	13/05/2022
18	ELSA2-AFP	29/04/2022	18/05/2022
18	ELSA-CA15-3	27/04/2022	18/05/2022
18	ELSA-CA19-9	27/04/2022	18/05/2022
18	ELSA-CA72-4	28/04/2022	18/05/2022
18	ELSA-FBHCG	28/04/2022	18/05/2022
17	MSK/TR	21/04/2022	18/05/2022
19	OCFE07-TRYPS	02/05/2022	18/05/2022
17	RBAN/25	21/04/2022	18/05/2022
19	THYRO	02/05/2022	18/05/2022
16	TR-AB-CT	11/04/2022	18/05/2022
19	TRYPS-DIA	02/05/2022	18/05/2022
19	TRYPS-US	02/05/2022	18/05/2022
16	REN-CT2P	15/04/2022	19/05/2022
19	BI-INS-IRMA	02/05/2022	20/05/2022
19	BI-INSU-US	02/05/2022	20/05/2022
18	ELSA-CYFRA-	28/04/2022	20/05/2022
19	ELSA-FBHCGSP	05/05/2022	20/05/2022
19	ELSA-NSE	03/05/2022	20/05/2022

SEMAINE de COMMANDE	REFERENCE KIT	DATE LIMITE de COMMANDE*	DEBUT des VENTES
19	OCFK07-PIIIP	04/05/2022	20/05/2022
19	PIIIP-US	04/05/2022	20/05/2022
19	AAI	03/05/2022	23/05/2022
19	AAI-US	03/05/2022	23/05/2022
19	CEA-RIACT	02/05/2022	23/05/2022
19	RENINE	03/05/2022	23/05/2022
18	ACHR-AB	27/04/2022	25/05/2022
19	ELSA-CA125II	04/05/2022	25/05/2022
20	ELSA-PTH	09/05/2022	25/05/2022
20	FPSA-RIACT	09/05/2022	25/05/2022
20	GASK-PR	09/05/2022	25/05/2022
20	GASK-PR-US	09/05/2022	25/05/2022
20	PSA-RIACT	09/05/2022	25/05/2022
18	TPO-AB-CT	27/04/2022	25/05/2022
20	CORT-CT2	12/05/2022	30/05/2022
20	CORT-US	12/05/2022	30/05/2022
20	ELSA-ACTH	11/05/2022	30/05/2022
20	ESTR-S-US	12/05/2022	30/05/2022
20	ESTR-US-CT	12/05/2022	30/05/2022
20	IRMA-HCT	12/05/2022	30/05/2022
20	OCFD03-FT4	11/05/2022	30/05/2022
20	OCFD03-FT4-K	11/05/2022	30/05/2022
20	OCFD53-FT4	11/05/2022	30/05/2022
20	OCFH07-FT3	11/05/2022	30/05/2022
20	OCFH57-FT3	11/05/2022	30/05/2022
20	OCPL07-TSH	10/05/2022	30/05/2022
20	OCPL57-TSH	10/05/2022	30/05/2022
19	TESTO-CT2	05/05/2022	30/05/2022
19	TESTO-US	05/05/2022	30/05/2022
19	TGAB I STEP	02/05/2022	02/06/2022
21	MSK/TR	19/05/2022	15/06/2022
21	RBAN/25	19/05/2022	15/06/2022
20	TR-AB-CT	12/05/2022	15/06/2022
21	17-OHP-CT	16/05/2022	16/06/2022
21	PROGEST-RIA	16/05/2022	16/06/2022
21	REN-CT2P	16/05/2022	16/06/2022
22	ACHR-AB	25/05/2022	22/06/2022
23	VGC/25	01/06/2022	29/06/2022
23	VGK/25	01/06/2022	29/06/2022
23	TGAB I STEP	30/05/2022	30/06/2022

* Au delà de ces dates de commande, merci de nous consulter pour connaître la disponibilité des produits correspondants.

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**

