Delegation granted for increase in capital by the Shareholders' meeting to the Board of Directors

Delegation of competence

Object - Duration	Limit	Use in 2020
Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right to the allotment of the debt instruments, with preferential subscription rights of the shareholders. (EGM 06/24/2020 – Resolution n°11)	The limit is $\pounds4,000,000$ corresponding to the maximum nominal amount of the increase of the share capital and to the maximal nominal amount of the debt instruments and $\pounds500,000,000$ on the maximum overall limit of the maximum nominal amount of the debt instruments.	None
Granted for a period of 26 months as from 24/06/2020	It being specified that the limits of the nominal amount of the capital increases and debt instrument, with or without preferential subscription rights of the shareholders, set from the twelfth (12 th) to the seventeenth (17 th) resolutions submitted to this Shareholders' Meeting shall be deducted from this overall limit	
Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right of the allotment of debt instruments, without preferential subscription rights of the shareholders – through public offerings,other than those referred to in the Article L. 411-2 of the French Monetary and Financial Code. (EGM 06/24/2020 – Resolution n° 12)	The limit is deducted on the overall limit of €4,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments).	None
Granted for a period of 26 months as from 24/06/2020		
Ability to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through public offers addressed exclusively to qualified investors or to a restricted circle of investors as defined in the article L. 411-2 of the French Monetary and Financial Code. (EGM 06/24/2020 – Resolution n° 13)	The limit is deducted on the overall limit of €4,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments).	None
Granted for a period of 26 months as from 24/06/2020		
Ability to increase the number of shares and/or securities giving access to the share capital of the Company to be issued in the event of a share capital increase with or without preferential subscription rights of the shareholders. (EGM 06/24/2020 – Resolution n° 14)	The limit amount 15% of initial issue of shares, pursuant to the resolution n°11 to 13 described above.	None
Granted for a period of 26 months as from 24/06/2020		
Ability to issue shares and/or securities giving access to the share capital of the Company, as consideration for securities tendered through public exchange offers initiated by the Company, without preferential subscription right of the shareholders.(EGM 06/24/2020 - Resolution n° 15)	share capital of the Company at the moment of the capital increase (increase of the share capital) and on the overall	None
Granted for a period of 26 months as from 24/06/2020		
Ability to increase the share capital through the capitalization of reserves, earnings or premiums or any other sum upon which capitalization would be permitted.	The limit is €4,000,000 (corresponding to the maximum nominal amount of the increase of the share capital); It is a independent limit.	None

Start 🔨

(EGM 06/24/2020 – Resolution n° 16)		
Granted for a period of 26 months as from 24/06/2020		
Ability to issue shares and/or securities giving access to the share capital giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders and reserved for members of saving plans. (EGM 06/24/2020 - Resolution n° 17)	The limit is €4,000,000 corresponding to the maximum nominal amount of the increase of the share capital; it is an independent limit.	None
Granted for a period of 26 months as from 24/06/2020		
Ability to reduce the capital by cancelling shares acquired under buyback program (FCM $0.6/24/2020$, Desclution $p^{\circ}(12)$)	The limit is of 10% of the capital of the Company and by period of 24 months.	None
(EGM 06/24/2020 - Resolution n°18)		
Granted for a period of 18 months as from 24/06/2020		
Ability to grant free new or existing shares to the benefit of employees or corporate officers	The limit amount of 10% of the Company's share capital calculated on the attribution date	None
(EGM 06/24/2020 - Resolution N°19)		
Granted for a period of 38 months as from 24/06/2020.		