

Aubagne, January 7, 2025

Half-year liquidity contract statement for Sartorius Stedim Biotech

Under the liquidity contract entered into between Sartorius Stedim Biotech and Kepler Cheuvreux, the following resources appeared on the liquidity account on December 31, 2024:

- 29,563 shares
- € 6,232,769.05
- Number of executions on buy side on semester: 5,450
- Number of executions on sell side on semester: 5,333
- Traded volume on buy side on semester: 281,035 shares for € 49,893,952.42
- Traded volume on sell side on semester: 297,764 shares for € 53,521,926.85

As a reminder :

The following resources appeared on the last half year statement on June 30, 2024 on the liquidity account:

- 46,292 shares
- € 2,521,189.91
- Number of executions on buy side on semester: 6,493
- Number of executions on sell side on semester: 5,322
- Traded volume on buy side on semester: 305,402 shares for € 69,412,550.00
- Traded volume on sell side on semester: 273,281 shares for € 63,664,027.43

The following resources appeared on the liquidity account when the activity started:

- 0 shares
- € 10,000,000.00

The implementation of this report is carried out in accordance with AMF Decision N°2021-01 of June 22, 2021 renewing the implementation of liquidity contracts for shares as an accepted market practice.

A profile of Sartorius Stedim Biotech

Sartorius Stedim Biotech is a leading international partner of the biopharmaceutical industry. As a provider of innovative solutions, the company based in Aubagne, France, helps its customers to manufacture biotech medications, such as cell and gene therapies, safely, rapidly, and economically. The shares of Sartorius Stedim Biotech S.A. are quoted on the Euronext Paris. The company has a strong global reach with manufacturing and R&D sites as well as sales entities in Europe, North America, and Asia. Sartorius Stedim Biotech regularly expands its portfolio through acquisitions of complementary technologies. In 2023, the company generated sales revenue of around 2.8 billion euros. Currently, more than 10,600 employees are working for customers around the globe.

Contact

Petra Müller

Head of Investor Relations

+49 (0)551.308.6035

Petra.mueller2@sartorius.com

Sartorius Stedim Biotech S.A.

Z.I. Les Paluds

Avenue de Jouques, CS 91051

13781 Aubagne Cedex, France

Phone: +33.(0)4.42.84.56.00

Fax: +33.(0)4.42.84.56.19

www.sartorius.com

	Buy side			Sell side		
	Number of executions	Number of shares	Traded volume in EUR	Number of executions	Number of shares	Traded volume in EUR
Total	5,450	281,035	49,893,952.42	5,333	297,764	53,521,926.85
7/1/2024	55	3,200	479,968.00	12	751	115,240.95
7/2/2024	79	4,448	641,312.64	84	4,500	658,080.00
7/3/2024	18	1,000	151,850.00	73	3,399	520,013.01
7/4/2024	19	1,400	217,784.00	62	2,700	422,739.00
7/5/2024	12	800	125,688.00	46	2,630	415,119.20
7/8/2024	41	2,300	365,999.00	33	1,700	272,017.00
7/9/2024	50	2,600	409,240.00	33	2,000	316,700.00
7/10/2024	20	1,100	173,305.00	30	1,700	270,045.00
7/11/2024	3	100	16,150.00	81	4,300	706,533.00
7/12/2024	39	2,017	336,496.11	25	1,841	311,975.86
7/15/2024	11	700	117,572.00	16	1,259	212,217.04
7/16/2024	11	500	83,760.00	37	3,000	506,610.00
7/17/2024	-	-	-	35	2,300	395,071.00
7/18/2024	-	-	-	38	3,100	549,661.00
7/19/2024	100	6,383	970,726.64	-	-	-
7/22/2024	45	2,900	421,544.00	6	300	44,349.00
7/23/2024	17	1,200	179,052.00	69	4,500	692,370.00
7/24/2024	61	3,400	532,780.00	59	3,800	603,136.00
7/25/2024	52	2,650	429,247.00	36	2,500	408,600.00
7/26/2024	7	500	82,120.00	77	3,900	664,482.00
7/29/2024	35	3,000	524,490.00	54	3,270	574,506.30
7/30/2024	65	4,300	748,114.00	106	6,800	1,201,220.00
7/31/2024	17	1,200	219,840.00	71	6,300	1,157,058.00
8/1/2024	31	2,100	387,114.00	50	3,000	556,320.00
8/2/2024	118	6,400	1,143,104.00	14	1,000	182,710.00
8/5/2024	80	5,000	853,900.00	34	2,400	412,920.00
8/6/2024	48	2,700	462,159.00	25	1,500	258,735.00
8/7/2024	14	1,500	262,545.00	40	3,100	544,732.00
8/8/2024	61	3,300	568,689.00	30	1,900	330,676.00
8/9/2024	35	1,700	299,149.00	31	1,600	283,888.00
8/12/2024	44	2,600	452,530.00	10	600	104,850.00
8/13/2024	53	3,600	602,388.00	46	2,600	442,390.00
8/14/2024	58	2,800	480,284.00	25	1,200	207,828.00
8/15/2024	30	1,600	271,568.00	33	2,000	341,780.00
8/16/2024	34	1,800	308,052.00	30	2,000	344,220.00
8/19/2024	46	1,300	224,445.00	32	1,970	341,873.80
8/20/2024	54	2,500	429,750.00	8	600	104,100.00
8/21/2024	4	400	68,696.00	40	2,100	362,040.00
8/22/2024	10	432	75,492.00	44	2,500	438,500.00
8/23/2024	35	1,900	331,189.00	40	2,200	386,496.00
8/26/2024	29	1,300	230,685.00	33	1,700	302,345.00
8/27/2024	19	1,000	179,070.00	48	2,700	485,298.00
8/28/2024	23	1,400	253,400.00	24	1,300	236,899.00
8/29/2024	20	963	174,794.13	32	2,200	402,512.00
8/30/2024	26	1,204	221,391.52	44	1,700	313,718.00
9/2/2024	47	2,683	477,252.04	-	-	-

	Buy side			Sell side		
	Number of executions	Number of shares	Traded volume in EUR	Number of executions	Number of shares	Traded volume in EUR
Total	5,450	281,035	49,893,952.42	5,333	297,764	53,521,926.85
9/3/2024	13	700	122,563.00	11	420	73,966.20
9/4/2024	50	2,108	358,360.00	25	1,400	241,276.00
9/5/2024	33	1,300	222,690.00	17	1,500	259,560.00
9/6/2024	28	1,900	339,872.00	53	3,500	628,285.00
9/9/2024	3	100	17,840.00	51	3,000	555,000.00
9/10/2024	18	1,420	270,183.40	28	1,600	307,216.00
9/11/2024	17	740	139,830.40	22	1,100	209,319.00
9/12/2024	40	1,480	279,320.40	26	1,000	190,880.00
9/13/2024	46	1,600	297,408.00	33	1,300	243,789.00
9/16/2024	69	2,800	511,028.00	34	1,800	330,048.00
9/17/2024	53	2,700	493,560.00	38	1,900	350,056.00
9/18/2024	80	4,600	810,796.00	71	3,900	709,371.00
9/19/2024	43	3,100	563,456.00	39	2,200	404,338.00
9/20/2024	86	4,823	841,468.81	-	-	-
9/23/2024	49	2,577	443,166.69	43	2,700	465,507.00
9/24/2024	34	2,300	396,727.00	31	1,850	320,327.50
9/25/2024	41	2,100	359,793.00	33	2,100	361,263.00
9/26/2024	-	-	-	90	5,900	1,060,525.00
9/27/2024	12	700	130,550.00	66	3,500	658,875.00
9/30/2024	49	3,000	573,390.00	37	2,322	447,426.18
10/1/2024	67	4,000	748,360.00	22	1,337	254,431.10
10/2/2024	55	2,800	512,232.00	18	1,000	183,900.00
10/3/2024	74	3,500	631,050.00	35	2,000	363,680.00
10/4/2024	46	2,370	428,116.80	42	2,400	439,272.00
10/7/2024	41	2,200	401,434.00	45	2,100	384,300.00
10/8/2024	55	2,514	448,472.46	45	2,600	467,636.00
10/9/2024	52	2,680	491,914.00	54	2,900	534,180.00
10/10/2024	46	2,400	437,688.00	55	2,800	513,324.00
10/11/2024	23	1,200	222,984.00	67	3,800	710,752.00
10/14/2024	24	1,600	302,000.00	52	3,083	586,139.96
10/15/2024	73	4,600	866,318.00	40	2,229	428,748.15
10/16/2024	92	3,886	690,464.48	55	2,500	451,100.00
10/17/2024	-	-	-	76	6,199	1,247,672.73
10/18/2024	2	100	20,800.00	63	3,900	821,886.00
10/21/2024	121	5,200	1,068,080.00	9	600	127,998.00
10/22/2024	101	4,300	852,819.00	72	3,734	755,238.84
10/23/2024	72	2,500	489,425.00	40	2,000	394,520.00
10/24/2024	65	3,500	678,930.00	26	1,600	312,352.00
10/25/2024	72	2,800	530,992.00	26	1,100	209,385.00
10/28/2024	43	2,471	469,761.81	64	2,600	496,340.00
10/29/2024	82	3,799	715,389.69	31	1,600	303,440.00
10/30/2024	86	3,500	643,475.00	46	2,700	497,799.00
10/31/2024	42	2,100	384,405.00	63	2,650	487,070.00
11/1/2024	12	600	109,548.00	49	2,800	524,664.00
11/4/2024	38	2,150	412,069.00	33	1,700	327,845.00
11/5/2024	27	1,550	292,128.50	21	1,100	208,098.00

	Buy side			Sell side		
	Number of executions	Number of shares	Traded volume in EUR	Number of executions	Number of shares	Traded volume in EUR
Total	5,450	281,035	49,893,952.42	5,333	297,764	53,521,926.85
11/6/2024	62	3,300	613,206.00	32	1,700	321,708.00
11/7/2024	10	700	129,311.00	22	1,400	259,756.00
11/8/2024	29	1,600	293,376.00	22	1,100	202,895.00
11/11/2024	25	1,200	220,764.00	37	1,400	258,146.00
11/12/2024	53	2,600	474,006.00	110	5,000	932,700.00
11/13/2024	64	3,400	622,200.00	43	2,000	368,600.00
11/14/2024	30	1,900	346,199.00	64	3,000	551,190.00
11/15/2024	143	6,500	1,139,970.00	6	400	70,760.00
11/18/2024	52	2,700	457,677.00	5	200	34,050.00
11/19/2024	73	3,300	548,097.00	44	2,117	354,682.18
11/20/2024	43	2,075	348,475.50	39	1,983	334,234.65
11/21/2024	57	2,390	396,644.40	54	2,700	449,928.00
11/22/2024	-	-	-	82	3,700	641,839.00
11/25/2024	27	1,700	301,223.00	55	3,300	587,070.00
11/26/2024	37	2,400	425,520.00	50	2,600	462,878.00
11/27/2024	34	2,000	359,920.00	56	2,200	397,188.00
11/28/2024	15	1,300	236,249.00	54	2,700	491,373.00
11/29/2024	38	1,950	350,883.00	32	1,800	324,558.00
12/2/2024	64	2,400	432,384.00	64	3,100	561,162.00
12/3/2024	61	2,900	514,344.00	21	1,000	177,670.00
12/4/2024	8	300	53,991.00	43	1,900	342,456.00
12/5/2024	75	3,825	683,030.25	16	1,100	199,298.00
12/6/2024	24	1,087	191,779.41	58	3,200	569,472.00
12/9/2024	38	1,600	288,176.00	68	3,400	617,100.00
12/10/2024	-	-	-	134	6,000	1,147,080.00
12/11/2024	46	1,717	328,204.55	51	3,400	653,004.00
12/12/2024	64	2,703	523,598.13	60	2,900	563,644.00
12/13/2024	46	2,200	427,900.00	40	2,100	410,340.00
12/16/2024	38	1,100	210,155.00	16	800	153,320.00
12/17/2024	70	2,940	558,012.00	33	2,200	420,640.00
12/18/2024	32	1,800	339,552.00	31	1,900	359,708.00
12/19/2024	90	3,801	706,986.00	38	2,200	411,378.00
12/20/2024	67	2,899	531,502.66	54	2,700	498,474.00
12/23/2024	19	1,000	184,780.00	20	1,300	240,916.00
12/24/2024	22	1,300	241,384.00	9	600	111,900.00
12/27/2024	12	800	149,360.00	45	2,520	472,147.20
12/30/2024	27	1,200	222,348.00	8	500	93,200.00
12/31/2024	4	200	37,200.00	24	1,200	224,088.00