

Sartorius Stedim Biotech S.A. - 13781 Aubagne, France

## Annual report of the liquidity contract with the stockbroker company Gilbert Dupont

Aubagne, France | January 3, 2020

This announcement contains corrections of the figures given in our annual report of the liquidity contract published yesterday.

Under the liquidity contract concluded between SARTORIUS STEDIM BIOTECH and the stockbroker Gilbert Dupont, the following assets appeared on the liquidity account on December 31, 2020:

- Number of shares: 3,225
- Liquidity account cash balance: €370,692.66

During the second half of 2019, the following were negotiated:

Buy	85,289 securities	€11,532,797.71	2,717 transactions
Sell	84,583 securities	€11,435,057.18	2,590 transactions

It should be noted that in the first half of 2019, which ended June 28, 2019, the following assets appeared on the liquidity account:

- Number of shares: 2,519
- Liquidity account cash balance: €468,433.16

It should also be noted that upon implementing the liquidity contract, the following assets were made available:

- Number of shares: 654
- Liquidity account cash balance: €394,895.12

### A profile of Sartorius Stedim Biotech

Sartorius Stedim Biotech is a leading international partner of the biopharmaceutical industry. As a total solutions provider, the company helps its customers to manufacture biotech medications safely, rapidly and economically. Headquartered in Aubagne, France, Sartorius Stedim Biotech is quoted on the Eurolist of Euronext Paris. With its own manufacturing and R&D sites in Europe, North America and Asia and an international network of sales companies, Sartorius Stedim Biotech has a global reach. The Group has been annually growing by double digits on average and has been regularly expanding its portfolio by acquisitions of complementary technologies. In 2018, the company earned sales revenue of €1,212.2 million and currently employs more than 6,000 people.

---

**Contact**

Ben Orzelek

Investor Relations

+49(0)551.308.1668

[ben.orzelek@sartorius.com](mailto:ben.orzelek@sartorius.com)

**Sartorius Stedim Biotech SA**

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-stedim.com](http://www.sartorius-stedim.com)

## ANNEXE S2 2019

	Buy			Sell		
	Number of trades	Number of shares traded	Turnover in EUR	Number of trades	Number of shares traded	Turnover in EUR
<b>TOTAL</b>	<b>2 717</b>	<b>85 289</b>	<b>11 532 797,71</b>	<b>2 590</b>	<b>84 583</b>	<b>11 435 057,18</b>
01/07/2019	16	600	82825,02	6	433	60443,29
02/07/2019	14	618	85323,98	12	585	81005,48
03/07/2019	6	306	42248,01	15	430	59726,01
04/07/2019	12	500	68765	0	0	0
05/07/2019	22	480	65417,52	6	165	22707,99
08/07/2019	7	300	40355,01	10	310	41994
09/07/2019	0	0	0	13	469	64029,38
10/07/2019	0	1020	139065,98	0	550	75104,98
11/07/2019	17	497	67704,22	20	797	109386,58
12/07/2019	20	684	93081,39	4	128	17574,4
15/07/2019	8	100	13660	18	1000	137790
16/07/2019	13	370	50832,01	34	1335	185016,05
17/07/2019	6	230	31980,99	10	762	106489,01
18/07/2019	9	316	43871,2	8	305	42511,51
19/07/2019	6	178	25899,39	12	276	40182,21
22/07/2019	4	300	44289,99	12	360	53228,99
23/07/2019	19	865	128936,04	25	950	142015,8
24/07/2019	25	781	116126,97	9	300	44990,01
25/07/2019	21	940	137290,02	1	30	4428
26/07/2019	6	200	29020	16	615	89674,5
29/07/2019	14	421	61206,92	6	210	30640,41
30/07/2019	11	386	55573,39	3	100	14532,5
31/07/2019	8	244	34883	10	365	52598,62
01/08/2019	10	255	36887,79	32	590	86070,09
02/08/2019	29	613	88490,41	14	298	43219,51
05/08/2019	47	997	137890,38	0	0	0
06/08/2019	14	259	34855,6	6	169	22849,59
07/08/2019	44	1499	202619,69	42	1084	146564,81
08/08/2019	16	420	56767,7	23	883	120608,09
09/08/2019	28	1088	147498,53	21	968	131451,6
12/08/2019	18	315	42595,21	15	328	44749,7
13/08/2019	11	295	39600,3	13	394	53371,52
14/08/2019	37	1667	224527,4	33	1467	197981,77
15/08/2019	18	376	49613,69	3	300	40074,99
16/08/2019	11	274	36393,69	37	962	128864,71
19/08/2019	2	28	3861,2	49	1082	147200,8
20/08/2019	24	516	71609,19	20	261	36493,59
21/08/2019	10	374	52096,4	29	511	71461,2
22/08/2019	46	1364	190902,7	2	50	7040
23/08/2019	19	379	52079	24	355	49112,69
26/08/2019	58	1470	198208,77	31	1462	197306,84
27/08/2019	37	955	128102,46	54	1707	229952,19
28/08/2019	28	586	80279,13	17	305	41911,09
29/08/2019	13	439	60028,42	24	520	71324,81
30/08/2019	20	514	71529,58	39	1168	162548,34
02/09/2019	27	804	112521,97	21	783	109976,73

03/09/2019	42	1137	158066,66	31	1126	156884
04/09/2019	1	25	3555	25	635	89668,41
05/09/2019	57	1488	208478,03	3	87	12421
06/09/2019	18	409	55985,88	3	90	12408
09/09/2019	37	667	89792,01	3	46	6248,8
10/09/2019	49	676	87928,87	23	429	55217,92
11/09/2019	9	319	41175,21	36	955	124401,83
12/09/2019	31	736	95309,72	32	941	122705,93
13/09/2019	15	521	67436,62	1	20	2612
16/09/2019	31	1359	173946,7	27	1031	132235,55
17/09/2019	4	150	19100	45	1458	188462,69
18/09/2019	11	440	57588,39	23	694	91051,13
19/09/2019	34	1240	163596,19	48	1677	221329,45
20/09/2019	58	1756	232699,84	22	809	107408,1
23/09/2019	38	910	119108,08	4	215	28159
24/09/2019	11	166	21358,51	35	1349	177151,62
25/09/2019	44	1279	164756,69	1	3	396
26/09/2019	16	153	19546,7	20	558	71962,08
27/09/2019	14	453	58197,82	18	632	81609,4
30/09/2019	27	905	116048,87	26	715	91880,07
01/10/2019	37	748	95340,6	22	623	79608,37
02/10/2019	27	600	74702,88	8	165	20560,01
03/10/2019	30	695	86413,38	30	781	97454,9
04/10/2019	17	709	88682,22	21	826	103725,2
07/10/2019	20	870	109077,9	26	888	111584,57
08/10/2019	31	1075	133796,66	12	920	114680,02
09/10/2019	33	1961	240319,38	30	2119	260153,88
10/10/2019	32	1474	180295,27	18	900	110210,04
11/10/2019	26	1555	188645,14	49	2276	277417,3
14/10/2019	9	250	30760	38	1708	211811,64
15/10/2019	11	311	39876,39	26	515	66298,01
16/10/2019	36	1111	143376,33	21	611	79256,17
17/10/2019	26	941	119269,49	13	662	84010,98
18/10/2019	32	722	91009,33	25	694	87496,33
21/10/2019	23	727	94582,41	66	1683	223865,6
22/10/2019	49	1961	258838,67	49	1420	188234,77
23/10/2019	50	965	124019,48	12	295	37495
24/10/2019	9	204	26104,8	39	1301	167875,19
25/10/2019	26	756	98866,81	17	722	94965,89
28/10/2019	17	562	73466,61	21	413	54417
29/10/2019	23	551	72161,99	10	430	56491,51
30/10/2019	10	285	37645,59	29	546	72706,02
31/10/2019	16	375	50216,59	12	311	41804,5
01/11/2019	18	809	108843,26	32	644	86799,99
04/11/2019	24	579	78791,07	28	521	71202,2
05/11/2019	21	651	88009,28	19	340	46054,5
06/11/2019	27	633	85186,23	25	648	87355,2
07/11/2019	7	461	62607,81	15	441	60397,42
08/11/2019	11	400	54160	11	365	49582,99
11/11/2019	12	584	79513,99	11	230	31475,59
12/11/2019	30	731	97701,51	15	358	47943,68
13/11/2019	5	92	12290,4	30	665	89818,03
14/11/2019	0	514	69519,48	0	511	69333,4

15/11/2019	18	519	70464,01	33	598	81441,98
18/11/2019	20	734	100296,92	22	780	107017,4
19/11/2019	18	582	79562,31	28	496	67947,98
20/11/2019	60	1839	249075,45	62	2119	287551,06
21/11/2019	33	1056	142230,95	10	347	46906,21
22/11/2019	9	372	49693,51	11	381	51314,99
25/11/2019	10	396	54913,72	42	873	120173,43
26/11/2019	3	250	35185	11	344	48981,99
27/11/2019	16	590	84350	23	629	90251,81
28/11/2019	34	867	123631,51	22	1159	165565,34
29/11/2019	6	300	42897,51	22	476	68313,9
02/12/2019	36	1091	155617,73	10	350	50767,5
03/12/2019	7	125	17537,5	25	1154	163315
04/12/2019	7	400	57170	17	632	90669,37
05/12/2019	13	815	118238,49	27	1178	171276,95
06/12/2019	19	665	96202,49	21	900	130802,49
09/12/2019	23	845	121944,49	6	450	64989,99
10/12/2019	33	1074	154100,42	16	658	94948,48
11/12/2019	17	417	59654,31	18	442	63500,42
12/12/2019	27	1325	190222,95	25	1143	164448,1
13/12/2019	49	1425	201884,88	10	454	65143,6
16/12/2019	0	0	0	8	267	38055
17/12/2019	15	411	59029,71	31	717	103220,68
18/12/2019	25	402	57855,4	6	159	23002,4
19/12/2019	4	50	7150	10	354	51029,7
20/12/2019	22	1150	167138,02	31	994	145045,08
23/12/2019	6	140	20538	11	529	77525,48
24/12/2019	6	250	37065	4	84	12398,2
27/12/2019	22	697	103066,71	6	256	37951,59
30/12/2019	13	343	50392,29	5	105	15486,5
31/12/2019	23	685	100606,43	12	431	63511,3

## ANNEXE S1 2019

	Number of trades			Number of shares traded		
	Number of trades	Number of shares traded	Turnover in EUR	Number of trades	Number of shares traded	Turnover in EUR
<b>TOTAL</b>	<b>2 404</b>	<b>72 168</b>	<b>7 778 802,43</b>	<b>2 522</b>	<b>73 065</b>	<b>7 889 953,10</b>
02/01/2019	59	2015	172365,92	43	1870	160639,73
03/01/2019	64	1676	141876,25	21	539	45994,7
04/01/2019	4	203	16956,35	49	1355	115196,41
07/01/2019	24	684	58504,5	29	822	70969,67
08/01/2019	19	617	53397,89	46	1152	100321,92
09/01/2019	29	926	82535,77	49	1134	101710,05
10/01/2019	31	720	64380,17	38	903	81450,33
11/01/2019	35	1225	110018,72	41	1039	93907,73
14/01/2019	56	1103	97355,52	0	0	0
15/01/2019	9	301	26386,41	31	739	65218,23

16/01/2019	20	564	49952,13	44	1349	120195,77
17/01/2019	46	1122	98755,64	0	0	0
18/01/2019	8	300	26352,51	27	1422	126139,93
21/01/2019	13	410	36250,07	27	881	78427,94
22/01/2019	30	1149	102148,86	20	539	48470,92
23/01/2019	30	1023	90858,77	11	471	42100,52
24/01/2019	13	525	46634,49	23	1057	94543,26
25/01/2019	3	100	8955	25	499	45184,45
28/01/2019	44	1313	117415,81	9	307	27565,74
29/01/2019	6	139	13781,35	75	2061	198471,62
30/01/2019	45	1550	154255,38	8	265	26725,01
31/01/2019	19	889	86310,52	8	200	19842,5
01/02/2019	21	786	76109,01	20	736	71754,11
04/02/2019	7	226	21906,5	15	498	48462,57
05/02/2019	6	200	19472,5	33	980	97098,5
06/02/2019	18	570	56723,32	16	745	74461,11
07/02/2019	18	428	42673,35	13	374	37452,51
08/02/2019	56	1072	105773,92	6	240	23810,5
11/02/2019	7	119	11793,55	29	864	86326,82
12/02/2019	48	1843	184321,38	36	936	94574,56
13/02/2019	0	0	0	43	1031	103711,49
14/02/2019	19	692	70449,61	25	579	59360,87
15/02/2019	12	700	70819	27	851	86898,93
18/02/2019	26	817	83280	30	782	79975,3
19/02/2019	19	791	80972,53	28	779	79956,4
20/02/2019	11	491	50335,5	36	799	82549,8
21/02/2019	22	936	97004,04	17	646	67391,62
22/02/2019	24	989	102400,76	31	1295	134205,64
25/02/2019	16	528	55247,91	23	806	84378,29
26/02/2019	28	1514	158305,05	19	1045	109495
27/02/2019	60	1372	142665,64	12	380	39697,5
28/02/2019	32	724	74731,5	26	581	60418,6
01/03/2019	0	0	0	42	952	99853
04/03/2019	0	0	0	13	378	40913,09
05/03/2019	14	411	44421,7	1	50	5430
06/03/2019	7	240	25776	24	486	52491,6
07/03/2019	19	513	55150,42	25	503	54351,72
08/03/2019	40	1052	112706,34	4	178	19108,41
11/03/2019	1	42	4452	14	434	46803,6
12/03/2019	5	94	10272,9	24	423	46358,39
13/03/2019	14	395	43929,02	18	571	63702,47
14/03/2019	28	714	80364,13	7	150	17050,01
15/03/2019	16	279	31288,4	18	272	30648,01
18/03/2019	10	399	44566,7	11	218	24409,7
19/03/2019	21	808	90726,52	23	743	83558,97
20/03/2019	34	770	86337,02	18	740	83171,12
21/03/2019	16	608	68113,51	16	464	52086,51
22/03/2019	22	675	75232,13	26	562	62818,28
25/03/2019	32	597	65335,38	17	561	61550
26/03/2019	2	200	21950	38	849	94565,02
27/03/2019	17	542	60566,82	7	102	11535,9
28/03/2019	28	622	69694,48	21	508	57169,81
29/03/2019	25	606	67573,73	41	998	111764,42

01/04/2019	12	171	19562,19	28	486	55472,38
02/04/2019	16	283	32553,6	10	180	20772,11
03/04/2019	3	49	5679,6	39	543	63044,09
04/04/2019	56	1312	151233,72	0	0	0
05/04/2019	21	1115	126046,18	23	782	88519,43
08/04/2019	11	475	53537,49	16	965	109234,91
09/04/2019	33	924	104238,57	20	504	57398,9
10/04/2019	16	348	38956,62	13	459	51668,12
11/04/2019	34	819	91727,59	42	1032	116024,46
12/04/2019	18	942	105436,74	23	975	109469
15/04/2019	20	725	81686,98	11	525	59296,28
16/04/2019	19	825	92680,01	22	600	67645,02
17/04/2019	22	796	87724,69	0	0	0
18/04/2019	0	0	0	32	1305	147261,16
19/04/2019	0	0	0	0	0	0
23/04/2019	6	355	37873,99	37	851	94121,96
24/04/2019	16	266	31016,61	34	748	87759,03
25/04/2019	22	654	78107,22	13	356	42663
26/04/2019	7	195	23205,6	7	235	28040,51
29/04/2019	14	606	72748,72	30	620	74475,21
30/04/2019	32	766	92555,47	33	853	103353,4
02/05/2019	25	554	66267,32	11	393	47318,89
03/05/2019	15	300	36080,01	13	477	57730,59
06/05/2019	13	523	62590,39	15	406	48863,6
07/05/2019	24	660	79657,12	21	692	84026,72
08/05/2019	12	337	40645,91	20	636	77389,12
09/05/2019	31	870	105720,31	8	324	39535,39
10/05/2019	10	373	45535,99	24	678	82853,7
13/05/2019	25	846	102312,19	2	15	1842
14/05/2019	6	155	18820,01	29	650	79644,11
15/05/2019	21	341	42299	17	422	52535,2
16/05/2019	8	225	27990	17	399	50293,91
17/05/2019	14	302	37961,1	12	304	38336,41
20/05/2019	18	516	64446,18	8	403	50784,21
21/05/2019	3	238	29617,01	7	244	30719,4
22/05/2019	11	615	77274,01	14	852	107394,17
23/05/2019	28	611	76994	5	416	52542,01
24/05/2019	23	1010	127100,02	2	200	25210
27/05/2019	7	215	27000	6	430	54180,99
28/05/2019	4	370	46220,99	13	750	94605
29/05/2019	9	614	76199,98	4	350	44019,99
30/05/2019	0	0	0	13	255	31530,5
31/05/2019	14	356	43768,81	16	575	70904,52
03/06/2019	14	507	62533,68	21	737	92809,6
04/06/2019	38	807	99460,57	23	637	78501,27
05/06/2019	3	164	20476,6	25	358	44887,9
06/06/2019	4	140	17738,99	8	179	22842,8
07/06/2019	5	141	18074,2	11	226	29180,19
10/06/2019	8	372	47831,61	8	164	21116,9
11/06/2019	7	149	19346,91	19	328	42807,12
12/06/2019	20	420	54664,89	10	330	43249,8
13/06/2019	14	389	50780,6	8	245	32075,99
14/06/2019	31	637	82148,6	8	257	33191,29

17/06/2019	15	430	55553,12	25	733	95385,51
18/06/2019	7	291	37978,79	25	440	58492,41
19/06/2019	14	500	67315,3	12	584	79094,51
20/06/2019	5	163	22356,49	16	219	30173,8
21/06/2019	42	858	118711,94	30	786	109447,2
24/06/2019	20	424	59112,38	6	105	14661,3
25/06/2019	22	501	69268,71	16	399	55526,6
26/06/2019	23	497	68005,01	8	365	50032,81
27/06/2019	5	202	27481,8	12	260	35621,61
28/06/2019	0	0	0	4	130	17978,51