



## Félicitations aux gagnants du 1<sup>er</sup> trimestre !

**Equipe L. Lejay,**  
INRA



**Equipe G. Bécard,**  
LRSV-CNRS 5546

**Equipe O. Loréal,**  
INSERM U991

CODE CHALLENGE HELIX	SORTIE JANV-14	SORTIE FEV-14	SORTIE MARS-14	TOTAL
12282	24	16	15	55
11577	14	12	9	35
11919	3	10	19	32
12276	12	11	6	29
12145	10	3	9	22
12420	3	7	11	21
12468	7	5	6	18
12392	3	3	7	13
12237	6	2	4	12
11567	1	6	5	12
11573	7	4	0	11
11745	5	2	3	10
11479	0	6	2	8
11585	3	3	2	8
12515	4	2	1	7
12097	1	2	4	7
11572	2	3	2	7
12525	2	2	2	6
12514	2	1	3	6
11923	2	2	2	6
12424	4	2	0	6
11967	1	3	1	5
11645	2	0	3	5
12320	3	1	1	5
12459	1	1	3	5
11653	1	1	1	3
12393	0	2	1	3
11801	0	3	0	3
12432	0	1	2	3
11642	1	0	1	2
11498	0	2	0	2
11652	0	0	1	1
11647	1	0	0	1
11749	0	1	0	1
12421	1	0	0	1
11664	0	0	0	0
12500	0	0	0	0
12314	0	0	0	0
12163	0	0	0	0