



Félicitations aux gagnants du 3^{ème} trimestre !

S. Borron,
INRA UMR BGP équipe EPI2V

Mr Loréal
INSERM U991
Hop. Pontchaillou



INSERM U1068

CHALLENGE HELIX	JUILLET 2014	Aout	Sept	Total Q3
11565	29	31	0	60
11919	19	2	13	34
12145	10	17	6	33
12276	9	1	5	15
12231	4	3	7	14
11573	10	1	2	13
11582	1	0	11	12
12514	4	4	3	11
12392	1	10	0	11
12421	9	2	0	11
11910	3	7	0	10
11745	3	1	5	9
12468	4	0	4	8
11923	6	1	1	8
12420	2	0	6	8
11585	1	4	3	8
11781	5	1	1	7
15480	7	0	0	7
12237	1	5	0	6
11967	1	1	4	6
12525	5	0	1	6
12441	1	1	4	6
11572	3	0	2	5
11652	1	2	2	5
12329	5	0	0	5
11839	3	1	1	5
11856	4	1	0	5
12097	1	3	0	4
11642	2	0	2	4
12459	3	0	1	4
12432	0	0	4	4
11554	0	1	3	4
12415	2	0	2	4
12435	1	1	2	4
12460	2	0	2	4
11922	2	1	1	4
11568	0	0	4	4

CODE CHALLENGE HELIX	SORTIE JUILLET 2014	Aout	Sept	Total Q3
12466	2	0	2	4
12433	1	2	1	4
11749	2	1	0	3
11567	1	0	2	3
12314	0	3	0	3
11838	2	0	1	3
12438	0	1	2	3
12440	1	1	1	3
12087	2	0	1	3
11479	2	0	0	2
11645	0	0	2	2
11829	0	0	2	2
11867	2	0	0	2
12416	0	0	2	2
11731	1	1	0	2
11882	1	0	1	2
12515	0	1	0	1
12320	1	0	0	1
11549	0	1	0	1
12431	0	0	1	1
11583	0	0	1	1
11653	0	0	0	0
11664	0	0	0	0
11647	0	0	0	0
12500	0	0	0	0
12393	0	0	0	0
11498	0	0	0	0
11801	0	0	0	0
12424	0	0	0	0
12163	0	0	0	0
12434	0	0	0	0
12404	0	0	0	0
12414	0	0	0	0
12409	0	0	0	0
12419	0	0	0	0

Le troisième Challenge est terminé. Votre ré-inscription est automatique à chaque nouveau Challenge. Toutefois, une seule chance pour remporter le bon Pass congrès est offerte par participant et pour l'année.