

Félicitations aux gagnants du 4^{ème} trimestre !

**C. BEAUDOIN, UMR GRED / PIAF,
Université Blaise Pascal,
Clermont Ferrand**

**Doan LUU,
INRA BPMP,
Montpellier**



**Mme BRIET, Laboratoire LBCMCP,
Université Toulouse
N. Prunier Leterme,
INRA Centre de Rennes, IGEPP, eq EGI, Le Rheu**



CODE CHALLENGE HELIX	SORTIE Oct 2014	Nov	Dec	Total
11972	16	13	11	40
12276	6	11	13	30
12420	11	10	2	23
12231	6	7	10	23
12083	10	6	6	22
12145	2	8	7	17
11919	7	2	3	12
11573	6	2	4	12
11583	3	4	4	11
12468	2	4	4	10
11554	4	1	4	9
11856	5	4	0	9
11582	4	1	4	9
11829	4	2	2	8
11479	4	1	2	7
12097	1	6	0	7
11745	2	2	3	7
12320	2	4	1	7
11585	2	2	3	7
11910	4	1	1	6
12460	4	2	0	6
12466	2	1	3	6
12316	4	0	2	6
12525	3	1	1	5
11923	1	1	3	5
12421	4	0	1	5
12329	2	2	1	5
11653	0	0	4	4
11749	1	0	3	4
11838	2	2	0	4
12438	1	3	0	4
11839	0	2	2	4
12087	2	2	0	4
12237	0	2	1	3
11572	0	2	1	3
11652	3	0	0	3

CODE CHALLENGE HELIX	SORTIE Oct 2014	Nov	Dec	Total
11567	2	1	0	3
12415	2	1	0	3
11922	0	0	3	3
11568	0	2	1	3
12515	1	1	0	2
11664	0	2	0	2
12514	0	1	1	2
11642	1	0	1	2
12459	2	0	0	2
12432	0	2	0	2
11549	0	0	2	2
11867	0	2	0	2
12431	0	2	0	2
12414	2	0	0	2
11967	0	0	1	1
12392	1	0	0	1
12314	0	0	1	1
12435	1	0	0	1
11781	0	1	0	1
12416	1	0	0	1
12404	0	0	1	1
11645	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
12440	0	0	0	0
12434	0	0	0	0
12441	0	0	0	0
15480	0	0	0	0
11731	0	0	0	0
12433	0	0	0	0
12409	0	0	0	0
11882	0	0	0	0
12419	0	0	0	0

1^{er} lot : 400€ pour un congrès scientifique

2^{ème} lot : Un bon d'achat de 100€ en produits Promega

3^{ème} lot : Un portoir timer & freezer

* Comptes remis à 0 pour le 3^{ème} Challenge. Réinscription automatique pour tous les participants du Challenge, exceptée pour l'équipe gagnante qui ne peut gagner plus d'une fois ces challenges.