

3^{ème} Challenge : Juillet 2014

RDV le **30 sept** pour connaître les gagnants

.... Et encore toutes nos félicitations à l'équipe de **M. Laouessergue** – gagnante du 2d Challenge



CODE CHALLENGE HELIX	SORTIE JUILLET 2014	Aout	Sept
11565	29		
11919	19		
12282	11		
12145	10		
11573	10		
12421	9		
12276	9		
11577	8		
15480	7		
11923	6		
12525	5		
12329	5		
11781	5		
12468	4		
12514	4		
11856	4		
11572	3		
11745	3		
12459	3		
11910	3		
11839	3		
11479	2		
11642	2		
11749	2		
12420	2		
12415	2		
11838	2		
12460	2		
11922	2		
11867	2		
12466	2		
12237	1		
11967	1		
12097	1		
12392	1		

CODE CHALLENGE HELIX	SORTIE JUILLET 2014	Aout	Sept
11652	1		
11567	1		
12320	1		
11585	1		
12435	1		
12440	1		
12441	1		
11731	1		
12433	1		
11582	1		
11653	0		
12515	0		
11664	0		
11645	0		
11647	0		
12500	0		
12393	0		
11498	0		
11801	0		
12314	0		
12424	0		
12432	0		
12163	0		
11549	0		
11554	0		
12438	0		
12434	0		
11829	0		
11568	0		
12416	0		
12404	0		
12431	0		
12414	0		
12409	0		
11583	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