

Penta C Allele Frequencies for Various Ethnic Groups

	African American		American Indian		Asian		Caucasian		Hispanic	
Allele	Number	Frequency	Number	Frequency	Number	Frequency	Number	Frequency	Number	Frequency
5	9	0.015	22	0.037	20	0.033	4	0.007	12	0.020
6	0	*	1	0.002	0	*	0	*	0	*
7	18	0.030	1	0.002	2	0.003	1	0.002	3	0.005
8	36	0.060	38	0.063	35	0.058	4	0.007	23	0.038
9	101	0.168	138	0.230	183	0.305	95	0.158	148	0.247
10	38	0.063	24	0.040	31	0.052	38	0.063	26	0.043
10.1	0	*	0	*	0	*	0	*	1	0.002
11	162	0.270	198	0.330	221	0.368	258	0.430	161	0.268
12	153	0.255	125	0.208	84	0.140	112	0.187	152	0.253
13	59	0.098	46	0.077	22	0.037	83	0.138	63	0.105
14	20	0.033	6	0.010	2	0.003	3	0.005	10	0.017
15	3	0.005	0	*	0	*	2	0.003	1	0.002
16	0	*	1	0.002	0	*	0	*	0	*
17	1	0.002	0	*	0	*	0	*	0	*
Sum	600	1.000	600	1.000	600	1.000	600	1.000	600	1.000
Persons Tested	300		300		300		300		300	
Heterozygosity	0.816		0.783		0.744		0.733		0.789	
HW Fischer	0.725		0.331		0.292		0.921		0.862	
PIC Value	0.790		0.751		0.705		0.696		0.756	
Matching Probability (pM)	0.022		0.032		0.047		0.054		0.029	
1/pM	44.450		31.161		21.407		18.470		34.411	
Probability of Discrimination	0.978		0.968		0.953		0.946		0.971	
Probability of Exclusion (A) (paternity)	0.639		0.583		0.525		0.515		0.587	
Typical Paternity Index (=1/(1-A))	2.766		2.397		2.106		2.063		2.421	
*Minimal allele frequency determinations are left to the user.										