

Y-STR Genotype.

Locus	Allele #1	Allele #2
DYS19	14	
DYS385a/b	13	16
DYS389I	14	
DYS389II	31	
DYS390	24	
DYS391	10	
DYS392	13	
DYS393	13	
DYS437	14	
DYS438	9	
DYS439	12	
DYS448	19	
DYS456	17	
DYS458	17	
DYS481	22	
DYS533	12	
DYS549	13	
DYS570	17	
DYS576	18	
DYS635	21	
DYS643	10	
Y-GATA-H4	11	