

**Population Genetics of 9 STRs in a Chinese  
Sample from Macau Using the  
“AmpF/STR™ Profiler Plus™ PCR Amplification Kit”**

Gusmão L<sup>1</sup>, SILVA F<sup>1</sup>, Alves C<sup>1</sup>, Amorim A<sup>1,2</sup>, Martins N<sup>3</sup>, Prata Mj<sup>1,2</sup>

<sup>1</sup> IPATIMUP

<sup>2</sup> Fac. Ciências, Universidade do Porto

<sup>3</sup> Centro de Transfusões de Sangue de Macau



Macau has been a Chinese territory under Portuguese administration for more than four centuries. A sample of unrelated individuals was typed for nine STR loci using “AmpF/STR™ Profiler Plus™ PCR Amplification Kit”.

Allele distributions were calculated for tetranucleotide markers: D3S1358, VW A, FGA, D8S1179, D21S11, D18S51, D5S818, D13S317 and D7S820. The data were analysed in terms of Hardy-Weinberg equilibrium and forensic informative value.

Estimated gene frequencies were:

D18S51	<b>13</b>	<b>14</b>	<b>15</b>	<b>16</b>	<b>17</b>	<b>18</b>	<b>19</b>	<b>20</b>	<b>21</b>	<b>22</b>	<b>23</b>	<b>24</b>
	12.0	21.3	21.3	17.6	7.4	5.6	4.6	3.7	2.7	1.9	0.9	0.9

D21S11	<b>28</b>	<b>29</b>	<b>29.2</b>	<b>30</b>	<b>30.2</b>	<b>31</b>	<b>31.2</b>	<b>32</b>	<b>32.2</b>	<b>33.2</b>
	1.9	29.6	0.0	25.9	1.9	8.3	12.0	2.8	12.0	5.6

FGA	<b>13</b>	<b>18</b>	<b>19</b>	<b>20</b>	<b>21</b>	<b>22</b>	<b>22.2</b>	<b>23</b>	<b>23.2</b>	<b>24</b>	<b>24.2</b>	<b>25</b>	<b>26</b>	<b>27</b>
	0.9	6.5	5.6	7.4	12.0	13.9	1.9	17.6	2.8	19.4	1.8	5.6	3.7	0.9

D8S1179	<b>8</b>	<b>10</b>	<b>11</b>	<b>12</b>	<b>13</b>	<b>14</b>	<b>15</b>	<b>16</b>	<b>17</b>
	0.9	16.7	12.0	8.4	17.6	18.5	14.8	10.2	0.9

vWA	<b>14</b>	<b>15</b>	<b>16</b>	<b>17</b>	<b>18</b>	<b>19</b>	<b>20</b>
	26.8	2.7	15.2	23.2	20.5	8.9	2.7

D7S820	<b>7</b>	<b>8</b>	<b>9</b>	<b>10</b>	<b>11</b>	<b>12</b>	<b>13</b>
	0.9	18.5	3.7	19.5	37.0	18.5	1.9