

Polymorphism of vWA locus in a family line

Toshinobu Komuro, Hirofumi Tsutsumi, Rei Mukoyama

Two types of structural polymorphism in vWA locus have been known, namely vWA-k (Kimpton et al.) and vWA-p (Ploos van Amstel et al.), respectively. In this report, we examined the vWA locus and found a new region which locates in between vWA-k and -p, and this locus was designated as vWA-t. The heredity relations between these three loci were also examined by using 23 DNA samples obtained from blood or oral mucosa (all of them are composed of four generations and belong to one family line). The DNA of the first generation was inherited to all other generations.