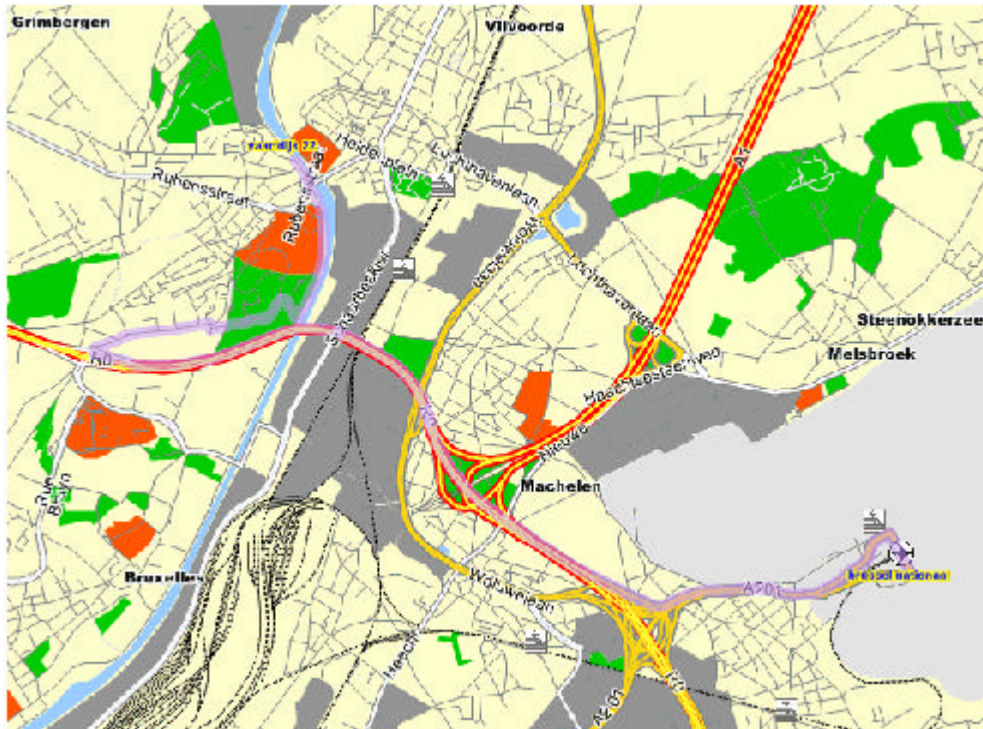
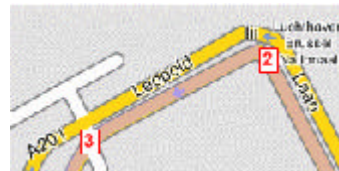


**FROM :** BRUSSEL NATIONAAL, ZAVENTEM, BELGIUM  
**TO :** VAARTDIJK 22, VILVOORDE, BELGIUM



- 1 Follow **Leopold III Laan / Vertrekhal** for 380m
- 2 Turn left, follow **Leopold III Laan** for 150m
- 3 At sign **UITRIT/SORTIE** turn left, follow the motorway **A201** for 1km
- 4 At sign **RING/(E411)/(E19)** exit : 4 Follow the motorway **R0-E40-E19** for 5km
- 5 At sign **KANAALZONE/MILITAIR ZIEKENHUIS** exit : 6 Follow **N216** for 450m



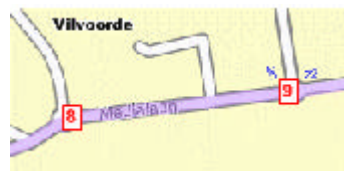
6 Turn right, follow **Koningslosteenweg** for 230m until **Vilvoorde**



7 Turn right, follow the roundabout for 30m



8 Turn right, follow **Medialaan** for 390m



9 Turn right, follow the roundabout for 30m



10 Turn right, follow **Beneluxlaan** for 260m



11 Turn right, follow **Drie Fonteinen** for 940m



12 Turn left, follow **N260** for 1km until **Vilvoorde**



13 Continue for 250m



14 Your reach **Vaardijk 22 Vilvoorde**



**FROM :** ANTWERP, BELGIUM  
**TO :** VAARTDIJK 22, VILVOORDE, BELGIUM

1 Follow the motorway **A1-E19** to **Brussels** for 30km

2 Follow the motorway **Vilvoorde** : exit **12** for 470m



3 Follow **N211 / Luchthavenlaan** for 1km



4 Turn right, follow the roundabout for 90m



5 Turn right, follow **N211** for 920m until **Vilvoorde**



- 6 Turn right, follow **Etienne Blondiastraat** for 150m
- 7 Turn left, follow **Van Helmontstraat** for 150m
- 8 Turn right, follow **N211 / Stationlei** for 250m
- 9 Follow **N1 / Heldenplein** for 790m
- 10 Follow **Rubensstraat** for 390m
- 11 Turn left, follow **N260 / Brusselsesteenweg** for 190m
- 12 Continue for 250m
- 13 Your reach **Vaardijk 22 Vilvoorde**

