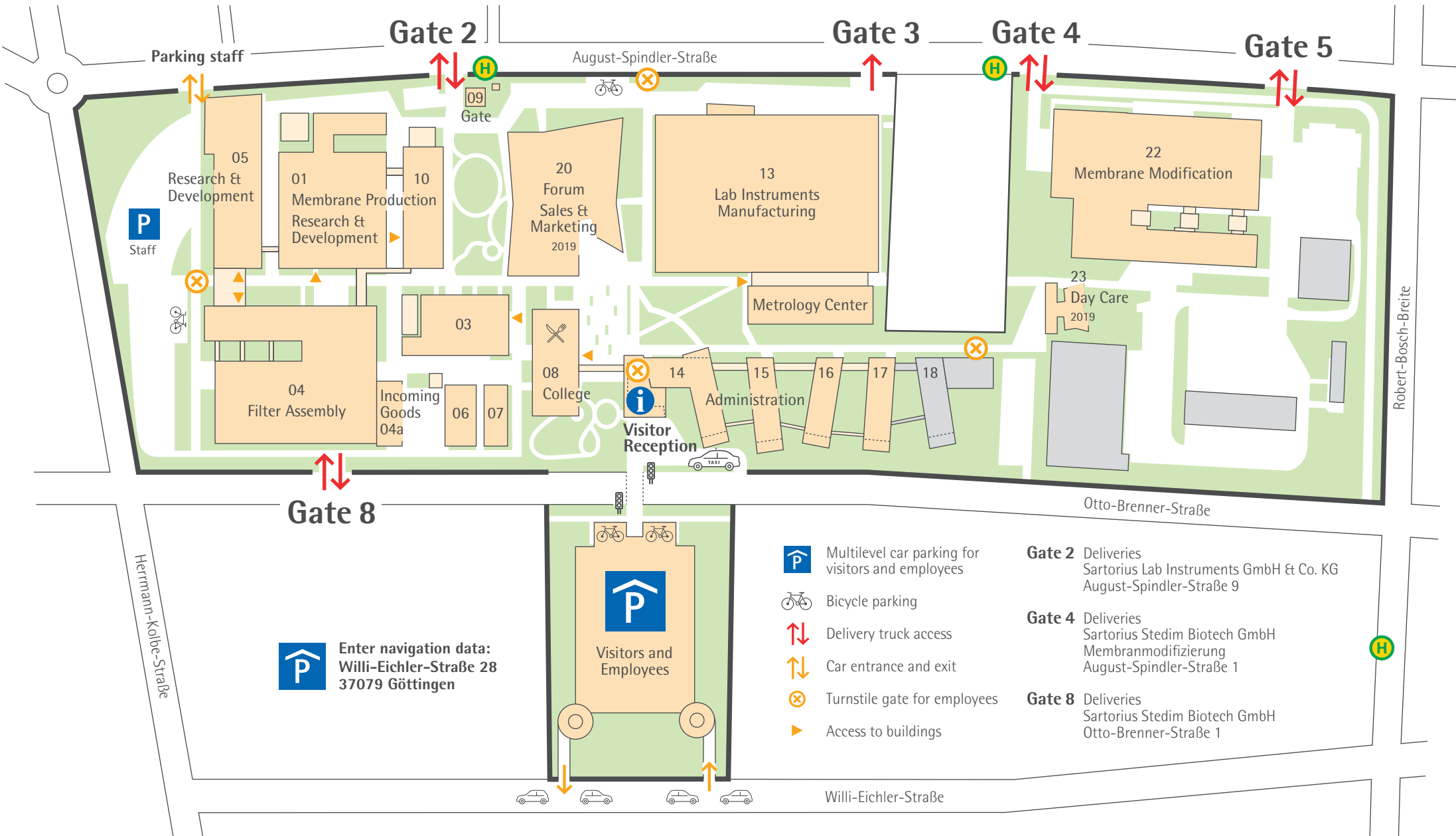



Map of Directions to Sartorius Campus



Access to Sartorius Campus



 Enter navigation data:
Willi-Eichler-Straße 28
37079 Göttingen

-  Multilevel car parking for visitors and employees
-  Bicycle parking
-  Delivery truck access
-  Car entrance and exit
-  Turnstile gate for employees
-  Access to buildings

- Gate 2** Deliveries
Sartorius Lab Instruments GmbH & Co. KG
August-Spindler-Straße 9
- Gate 4** Deliveries
Sartorius Stedim Biotech GmbH
Membranmodifizierung
August-Spindler-Straße 1
- Gate 8** Deliveries
Sartorius Stedim Biotech GmbH
Otto-Brenner-Straße 1