

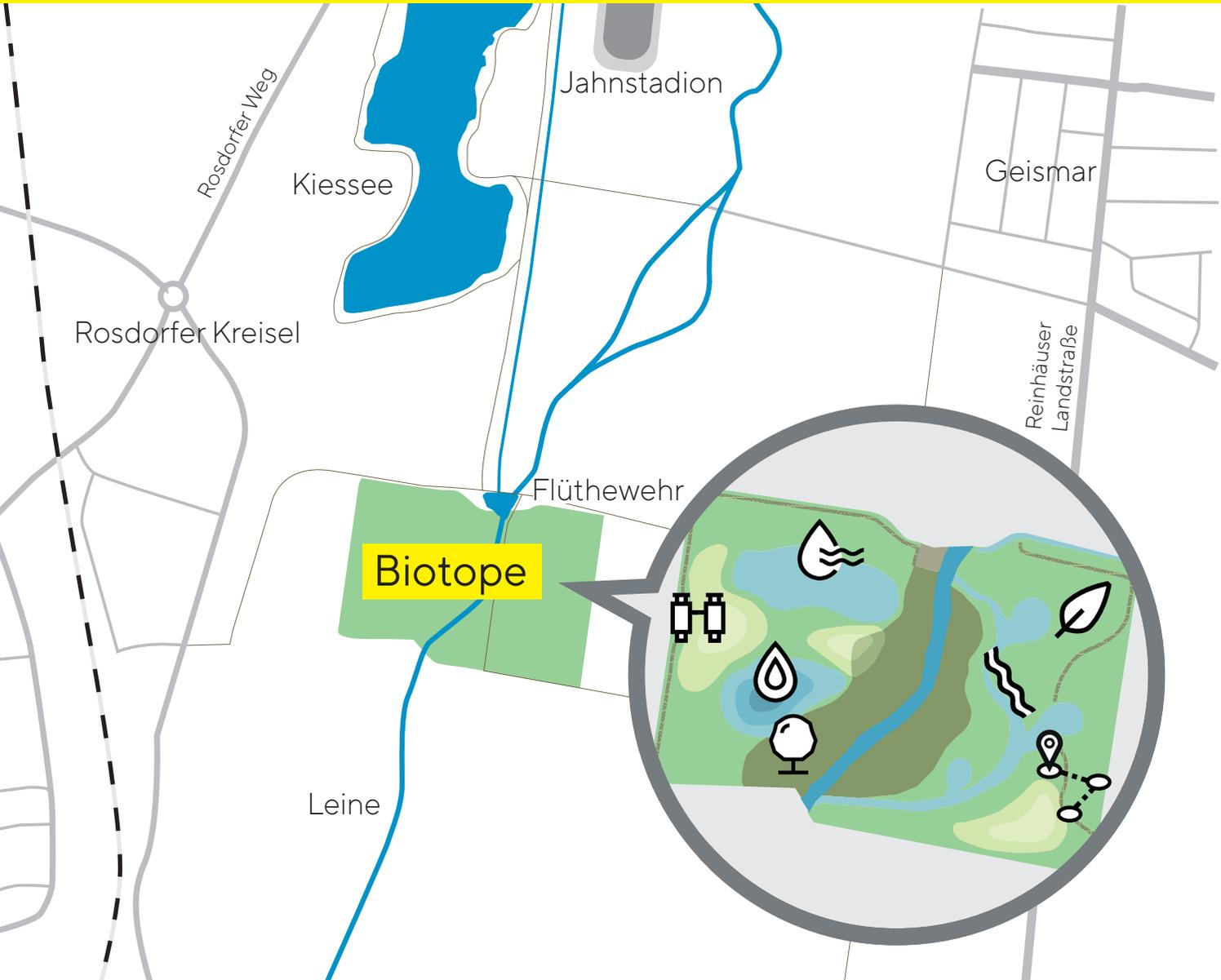
Biotope at the Kieselsee

Sartorius launched an initiative to create a diverse wetland biotope south of Flüthwehr near the Kieselsee recreational area in Göttingen by autumn 2021. The Heinz Sielmann foundation will coordinate the project, and the city of Göttingen will provide a 16-hectare area – around 40 acres – at the Leine River, previously used for predominantly agricultural purposes. Part of this area is on city district territory. Leine River divides the area into two even sections, which will provide an ecologically diverse habitat for rare and endangered animals, such as amphibians and birds.

Start of preparations:
February 2021



Planned opening:
Autumn 2021



-  Shallow pond
-  Observation point
-  Pond
-  Riparian forest
-  Extensive grassland
-  Flood channel
-  Experience trail

Heinz
Sielmann
Stiftung

SARTORIUS