



Feuchtbiotop

-  Flachwassertümpel
-  Beobachtungsturm
-  Stillgewässer
-  Weichholz-Auwald
-  Extensivgrünland
-  Flutrinne mit Mulden
-  Erlebnisweg

Heinz
Sielmann
Stiftung

SARTORIUS