

# How to find us



## By car:



**Frankfurt Airport (Germany) to Bad Wildbad:** approx. 2.5 hours  
 Distance: Approx. 180 km  
 Frankfurt Airport: Autobahn **A5** to Karlsruhe  
 Karlsruhe: Change to Autobahn **A8** direction Pforzheim/Stuttgart  
 Pforzheim: Take exit Pforzheim West, Exit 43!  
 B294: Follow the federal road **B294** to Birkenfeld → Neuenbürg → Höfen → Calmbach  
 At Calmbach: Leave **B294** at the entrance to Calmbach and follow the turn-off to → Bad Wildbad/Enzklösterle

**Stuttgart Airport (Germany) to Bad Wildbad:** approx. 1.5 hours  
 Distance: Approx. 70 km  
 Stuttgart airport: Autobahn **A8** to Pforzheim  
 Pforzheim: Take exit Pforzheim West, Exit 43!  
 B294: Follow the federal road **B294** to Birkenfeld → Neuenbürg → Höfen → Calmbach  
 At Calmbach: Leave **B294** at the entrance to Calmbach and follow the turn-off to → Bad Wildbad/Enzklösterle

**Strasbourg Airport (France) to Bad Wildbad:** approx. 1.5 hours  
 Distance: Approx. 120 km  
 Strasbourg airport: Federal road direction Offenburg take Autobahn **A5** direction Karlsruhe  
 Karlsruhe: Autobahn **A8** direction Pforzheim/Stuttgart  
 Pforzheim: Take exit Pforzheim West, Exit 43!  
 B294: Follow the federal road **B294** to Birkenfeld → Neuenbürg → Höfen → Calmbach  
 At Calmbach: Leave **B294** at the entrance to Calmbach and follow the turn-off to → Bad Wildbad/Enzklösterle

## By train:



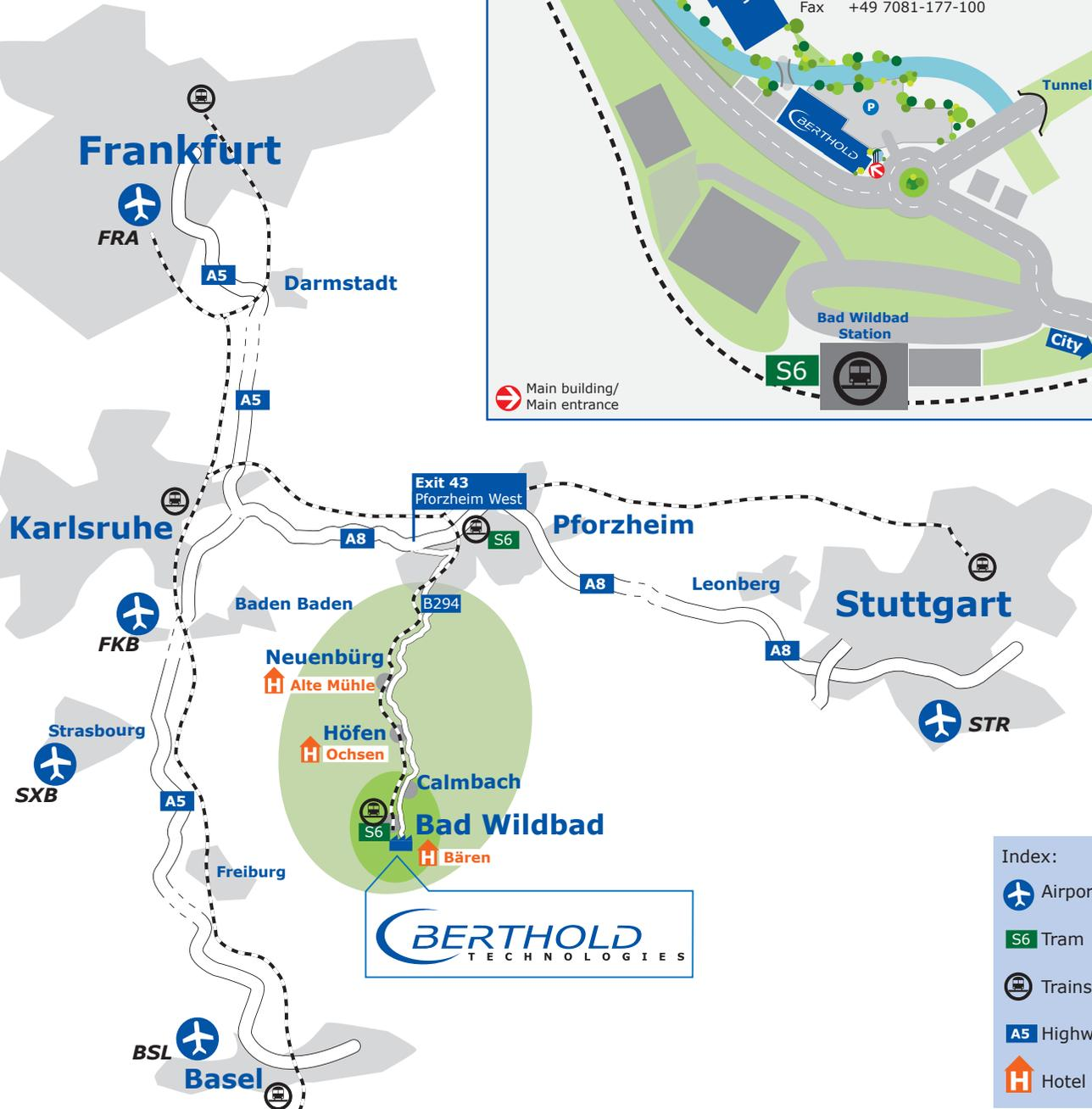
**Frankfurt Airport (Germany) to Bad Wildbad:** approx. 2.5 hours  
 Frankfurt airport station to Karlsruhe Railway station  
 Main station Karlsruhe to Pforzheim  
 Main station Pforzheim Tram **S6** to Bad Wildbad

**Stuttgart Airport (Germany) to Bad Wildbad:** approx. 2 hours  
 Subway station Stuttgart airport to Stuttgart Main station  
 Main station Stuttgart to Pforzheim  
 Main station Pforzheim Tram **S6** to Bad Wildbad

## Coordinates:



48° 45' 23.30" N  
 8° 33' 09.19" O  
 Altitude: 422 m NN



### Index:

- Airport
- S6 Tram
- Trainstation
- A5 Highway
- Hotel