

Selection of Hotels in Bad Wildbad

| Name and Address | Phone/Fax/eMail/Internet | 2019 Prices in € per person/night S – single room D- double room |
|---|--|--|
| Mokni's Palais Hotels & Spa Hotel Rossini und Badhotel Am Kurplatz 4-6 75323 Bad Wildbad | Phone +49 7081 301-0 Fax +49 7081 301-166 eMail: hotel@moknis.com Internet: www.moknis.com Prices include breakfast buffet and entrance to Palais Thermal. American Express card is not accepted | Rossini S 86,00 Rossini D 82,00 Badhotel S 105,00 Badhotel D 100,00 |
| Hotel Weingärtner Olgastraße 15 75323 Bad Wildbad | Phone +49 7081 1706-0 Fax +49 7081 1706-70 eMail: weingaertner-hotel@t-online.de Internet : www.hotel-weingaertner.com Prices include breakfast buffet and use of wellness area | S 65,00 D 57,00 |
| Hotel Rothfuß Olgastraße 47 75323 Bad Wildbad | Phone +49 7081 92 48 0 Fax +49 7081 92 48 10 eMail: anfrage@hotel-rothfuss.de Internet: www.wellnesshotel-rothfuss.de Prices include breakfast buffet and use of wellness area | S 78,00 S 78,00 |
| Hotel Auerhahn Heermannsweg 4 75323 Bad Wildbad (Sommerberg) | Phone +49 7081 95213 50 eMail: info@auerhahn-badwildbad.de Internet: www.auerhahn-badwildbad.de Prices include breakfast buffet | S 59,00 D 82,00 |
| Hotel Bergfrieden Bätznerstr. 78 75323 Bad Wildbad | Phone +49 7081 1704-0 Fax +49 7081 1704-20 eMail: empfang@hotelbergfrieden.de Internet: www.hotelbergfrieden.de Prices for Comfort plus room incl. breakfast | S 55,10 D 48,45 |
| Gasthof Zur Alten Linde Wilhelmstraße 74 75323 Bad Wildbad | Phone +49 7081 926-0 Fax +49 7081 926-250 Prices include breakfast buffet | S 50,00 D 50,00 |

**Please state that you are a visitor of Berthold Technologies when booking the room
 and you will receive a special discount.**

| Name and Address | Phone/Fax/eMail/Internet | 2019 Prices in € per person/night S – single room D- double room |
|---|---|---|
| Hotel Sonnenhof Bismarckstrasse 23 75323 Bad Wildbad | Phone +49 7081 92960 Fax +49 7081 929696 eMail: info@sonnenhof-badwildbad.de Internet: www.sonnenhof-badwildbad.de Prices incl. breakfast buffet and use of wellness area | S 59,00 D 53,00 |

Selection of Hotels in the region

| Name and Address | Phone/Fax/eMail/Internet | 2019 Prices in € per person/night S – single room D- double room |
|--|--|---|
| Hotel Ochsen Bahnhofstraße 2 75339 Höfen (reachable per car & tram – about 6 kms to Bad Wildbad) | Phone +49 7081 791-0 Fax +49 7081 791-100 eMail: info@ochsen-hoefen.de Internet: www.ochsen-hoefen.de Prices for business rooms incl. breakfast buffet and use of wellness area | S 70,05 D 54,75 |
| Hotel zur Alten Mühle Im Gänzbrunnen 75305 Neuenbürg (reachable per car and tram- about 8 kms to Bad Wildbad) | Phone +49 7082 924-00 Fax +49 7082 924-099 eMail: zordel@t-online.de Internet: www.zordel.de Prices include breakfast buffet (free WLAN) | S 79,00 D 60,50 |
| Berlin's Krone Lamm Marktplatz 2+3 75385 Bad Teinach-Zavelstein (reachable per car – about 17 kms from Bad Wildbad) | Phone +49 7053 9294-0 Fax +49 7053 9294-30 eMail: kronelamm@berlins-hotel.de Internet: www.berlins-hotel.de Prices include breakfast buffet and use of wellness area | Krone: S 95,00 Lamm: S 129,00 |

**Please state that you are a visitor of Berthold Technologies when booking the room
and you will receive a special discount.**