

Selection of Hotels in Bad Wildbad

Name and Address	Phone/Fax/eMail/Internet	2018 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. Credit cards are only accepted from Berthold customers. No Wi-Fi in the room.	Rossini S 85,00 Rossini D 81,00 Badhotel S 92,00 Badhotel D 88,00
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 63,00 D 55,00
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 79,00 from 3 nights S 75,00
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 include breakfast and free entrance into the Vital Therme	S from 41,80 D from 35,15
Gasthof Zur Alten Linde Wilhelmstraße 74 75323 Bad Wildbad	Phone +49 7081 926-0 Fax +49 7081 926-250 Prices include breakfast buffet	S 53,00 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	2018 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 include breakfast buffet and use of wellness area	S 55,00 D 52,00

Selection of Hotels in the region

Name and Address	Phone/Fax/eMail/Internet	2018 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 incl. breakfast buffet and use of indoor pool / sauna (DSL connection with cable free of charge, payable via hotspot)	S from 70,05 D from 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, use of wellness area, 0.5 l mineral water p. P./day in the room, tourist tax	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.**