

# Nanion's 9<sup>th</sup> Annual User Meeting

## When:

October 2018  
Start: Thursday 25<sup>th</sup> 1:00 PM  
End: Friday 26<sup>th</sup> 3:00 PM

## Where:

Nanion headquarters  
Ganghoferstraße 70a  
80339 Munich

## Arriving at Munich main station:

Take subway U5 (every 10 min)  
get off at Schwanthalershöhe  
and walk to Ganghoferstr. 70a (400 m).  
Duration approx. 10 min.



## Registration:

The User Meeting is free to attend. Nanion does not cover any travel or hotel expenses.



## Presentation:

If you like to present your data during the poster session or symposium: Please write an e-mail to: [gesa.@nanion.de](mailto:gesa.@nanion.de)

## Arriving at Munich airport:

Take train S8 (every 20 min), get off at Donnersberger Brücke. From here, take the Bus #53 direction Aidenbachstraße and get off at Ridlerstraße. Walk to Ganghoferstr. 70a (80 m). Duration: approx. 1h

# Nanion's 9<sup>th</sup> Annual User Meeting

## Hotels:

### 1. Sheraton Munich Westpark Hotel

Garmischer Str. 2, 80339 Munich

<http://www.sheratonwestpark.com/en>

10 minutes walking distance to headquarters  
- Single room: 129,00 Euro including breakfast

25 rooms are reserved for you until September 13.  
Please call +49 89 5196-6750 and book your room  
Mention: „ Nanion User Meeting 2018“

User Meeting Dinner will take place at the Seraton Hotel Restaurant



### 2. Hotel Ars Vivendi

Anglerstr. 19, Schwanthaler Höhe, 80339 Munich

<http://www.hotel-ars-vivendi.de/>

5 minutes walking distance to headquarters

All prices include breakfast:

- Double room standard: 1 person 110,00 Euro

12 Rooms are reserved for you until September 13.  
Please write an e-mail to: [info@hotel-ars-vivendi.de](mailto:info@hotel-ars-vivendi.de)  
Enter in reference line: „ Nanion User Meeting 2018“



*Ars Vivendi*  
\*\*\* HOTEL \*\*\*

### 3. Hotel Rivoli

Albert-Rosshaupter-Str. 18, 81369 Munich

<https://hotelrivoli.de>

10 minutes walking/ subway distance to  
headquarters

- Single room: 95,00 Euro including breakfast

15 rooms are reserved for you until September 13.  
Please write an e-mail to: [info@rivoli.de](mailto:info@rivoli.de)  
Enter in Reference line: „ Nanion User Meeting 2018“



HOTEL RIVOLI  
★★★ S  
MÜNCHEN

