

Nanion's 8th Annual User Meeting

When:

October 2017

Start: Wednesday 25th 1:00 PM

End: Thursday 26th 2:00 PM

Where:

Nanion headquarters
Ganghoferstraße 70a
80339 Munich



Arriving from Munich main station:

Take subway U5 (every 10 min)
get off at Schwanthalerhöhe 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.



Arriving from 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 (77m).
Duration: approx. 1h

Presentation:

If you like to present your data during the
poster session or symposium:
Please write an e-mail to:
gesa.rascher-eggstein@nanion.de

Nanion's 8th Annual User Meeting

Hotels:

1. 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:

- Single room standard: 82,00 Euro
- Single room comfort: 88,00 Euro
- Double room standard: 1 person: 93,00 Euro; 2 persons 127,00 Euro
- Double room comfort: 1 person: 98,00 Euro; 2 persons 132,00 Euro

28 Rooms are reserved for you.

Please write an e-mail to: Rico.Stiehl@hotel-ars-vivendi.de

Enter in reference line: „ [Nanion User Meeting 2017](#)“



2. Bavaria Boutique Hotel

Gollierstr. 9, 80339 Munich

<https://www.hotel-bavaria.com/de/>

15 minutes walking distance to headquarters

All prices include breakfast:

- Single room: 123,00 Euro

15 Rooms are reserved for you.

Please write an e-mail to: info@hotel-bavaria.com

Enter in reference line: „ [Nanion User Meeting 2017](#)“

**BAVARIA
BOUTIQUE HOTEL**



3. Sheraton Munich Westpark Hotel

Garmischer Str. 2, 80339 Munich

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

10 minutes walking distance to headquarters

Please book directly via Sheraton's homepage.



Nanion's 8th Annual User Meeting

Maps:



If you have questions, please do not hesitate to send an e-mail to:
gesa.rascher-eggstein@nanion.de
I am happy to help.

We are looking forward to seeing you soon!