

Nanion's Grand Copenhagen Opening Party

Nanion is proud to announce the opening of its Copenhagen Company Site





Friends, customers, neighbors, partners, Danes and curious are highly welcome to celebrate with us!

Monday, July 4th 18:00 o'clock Café Væksthuset at Frederiksberg Campus, Grønnegårdsvej 1 (55°40'51.3"N 12°32'25.9"E)

FENS ATTENDEES:

A bus shuttle will pick you up at 18:15 o'clock from The Bella Center Front entrance

Don't miss:

- Nanion's Lab tour
- Greenhouse Dinner