

Impressions from the Grand Opening
Celebration of our New Company Site in
Copenhagen (4th of July 2016)



„We would like to thank our clients and friends who joined our celebration and made this evening a wonderful and unforgettable evening for us. Your Nanion team.“

Impressions from the Grand Opening
Celebration of our New Company Site in
Copenhagen (4th of July 2016)



„We would like to thank our clients and friends who joined our celebration and made this evening a wonderful and unforgettable evening for us. Your Nanjion team.“