



complexity simplified.

## Luminex Corporation to Present at 39th Annual J.P. Morgan Healthcare Conference

January 13, 2021

AUSTIN, Texas, Jan. 13, 2021 /PRNewswire/ -- As previously announced, Luminex Corporation (Nasdaq: LMNX) will present at the 39<sup>th</sup> Annual J.P. Morgan Healthcare Conference, to be held virtually today, January 13, 2021 at 3:30 p.m. Central Time. A live webcast of the Luminex presentation has been made available and may be accessed at the company's website at <https://investor.luminexcorp.com/events-presentations>. The presentation will be available on our website within 24 hours and archived for six months on the website using the 'replay' link.

### About Luminex Corporation

At Luminex, our mission is to empower labs to obtain reliable, timely, and actionable answers, ultimately advancing health. We offer a wide range of solutions applicable in diverse markets including clinical diagnostics, pharmaceutical drug discovery, biomedical research, genomic and proteomic research, biodefense research, and food safety. We accelerate reliable answers while simplifying complexity and deliver certainty with a seamless experience. To learn more about Luminex, please visit us at [www.luminexcorp.com](http://www.luminexcorp.com).

#### Contact:

Harriss Currie  
Senior Vice President and Chief Financial Officer  
512-219-8020  
[hcurrie@luminexcorp.com](mailto:hcurrie@luminexcorp.com)

Carla Stanaford  
Investor Relations  
937-469-2120  
[cstanaford@luminexcorp.com](mailto:cstanaford@luminexcorp.com)



 View original content to download multimedia:<http://www.prnewswire.com/news-releases/luminex-corporation-to-present-at-39th-annual-jp-morgan-healthcare-conference-301207679.html>

SOURCE Luminex Corporation