



complexity simplified.

Luminex Corporation To Release Third Quarter 2019 Results After Market Close On November 4, 2019

October 25, 2019

AUSTIN, Texas, Oct. 25, 2019 /PRNewswire/ -- Luminex Corporation (NASDAQ: LMNX) today announced that it plans to report results for the third quarter ended September 30, 2019 on Monday, November 4, 2019 after the market close.

The company will host a conference call that day at 5:00 p.m. Eastern Time to discuss operating highlights and financial results.

The conference call will be webcast live and may be accessed at Luminex's Investor Relations website at <https://investor.luminexcorp.com>. The presentation slides will be posted to our Investor Relations website after the market close on November 4, 2019. Analysts and institutional investors may participate on the conference call by dialing (877) 930-7053 (U.S.) or (253) 336-7290 (outside the U.S.). The access code is 1293259. The webcast will be archived for six months on our 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.

Contacts:

Luminex Corporation

Jeff Christensen

Senior Director, Investor Relations

jeffc@luminexcorp.com

512-249-3033

Lazar FINN Partners

David Carey

david.carey@finnpartners.com

212-867-1768



 View original content to download multimedia: <http://www.prnewswire.com/news-releases/luminex-corporation-to-release-third-quarter-2019-results-after-market-close-on-november-4-2019-300945751.html>

SOURCE Luminex Corporation