



complexity simplified.

## Luminex Corporation Declares Third Quarter Cash Dividend

September 15, 2020

AUSTIN, Texas, Sept. 15, 2020 /PRNewswire/ -- On September 10, 2020, Luminex Corporation (Nasdaq:LMNX) (the "Company"), announced that its board of directors declared a cash dividend for the third quarter of 2020 of \$0.09 per share of common stock payable on October 15, 2020 to stockholders of record as of the close of business September 24, 2020.

### 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-declares-third-quarter-cash-dividend-301130669.html>

SOURCE Luminex Corporation