



complexity simplified.

## Luminex Corporation Declares First Quarter Cash Dividend

February 8, 2021

AUSTIN, Texas, Feb. 8, 2021 /PRNewswire/ -- Luminex Corporation (Nasdaq:LMNX) (the "Company") today announced that its board of directors declared a cash dividend for the first quarter of 2021 of \$0.10 per share of common stock payable on April 15, 2021 to stockholders of record as of the close of business March 25, 2021.

### 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).

### Contacts:

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

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-first-quarter-cash-dividend-301224092.html>

SOURCE Luminex Corporation