



complexity simplified.

Luminex Corporation Declares Second Quarter Cash Dividend

May 20, 2021

AUSTIN, Texas, May 20, 2021 /PRNewswire/ -- Luminex Corporation (Nasdaq:LMNX) (the "Company") today announced that its board of directors declared a cash dividend for the second quarter of 2021 of \$0.10 per share of common stock payable on July 8, 2021 to stockholders of record as of the close of business June 17, 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.

Contacts:

Harriss Currie
Senior Vice President of Finance and
Chief Financial Officer
hcurrie@luminexcorp.com
512-219-8020

Carla Stanaford
Investor Relations
937-469-2120
cstanaford@luminexcorp.com

 View original content to download multimedia: <http://www.prnewswire.com/news-releases/luminex-corporation-declares-second-quarter-cash-dividend-301296427.html>

SOURCE Luminex Corporation