

UNITED STATES
SECURITIES AND EXCHANGE COMMISSION
Washington, D.C. 20549

SCHEDULE 14A

Proxy Statement Pursuant to Section 14(a) of the
Securities Exchange Act of 1934 (Amendment No.)

Filed by the Registrant

Filed by a Party other than the Registrant

Check the appropriate box:

- Preliminary Proxy Statement
- Confidential, for Use of the Commission Only (as permitted by Rule 14a-6(e)(2))**
- Definitive Proxy Statement
- Definitive Additional Materials
- Soliciting Material Pursuant to §240.14a-12

LUMINEX CORPORATION

(Name of Registrant as Specified In Its Charter)

(Name of Person(s) Filing Proxy Statement, if other than the Registrant)

Payment of Filing Fee (Check the appropriate box):

- No fee required.
- Fee computed on table below per Exchange Act Rules 14a-6(i)(1) and 0-11.

- (1) Title of each class of securities to which transaction applies:
- (2) Aggregate number of securities to which transaction applies:
- (3) Per unit price or other underlying value of transaction computed pursuant to Exchange Act Rule 0-11 (set forth the amount on which the filing fee is calculated and state how it was determined):
- (4) Proposed maximum aggregate value of transaction:
- (5) Total fee paid:

- Fee paid previously with preliminary materials.
- Check box if any part of the fee is offset as provided by Exchange Act Rule 0-11(a)(2) and identify the filing for which the offsetting fee was paid previously. Identify the previous filing by registration statement number, or the Form or Schedule and the date of its filing.

- (1) Amount Previously Paid:
- (2) Form, Schedule or Registration Statement No.:
- (3) Filing Party:
- (4) Date Filed:



LUMINEX CORPORATION
ATTN: KENDEL MARTIN
12212 TECHNOLOGY BLVD.
AUSTIN, TX 78727

Your **Vote** Counts!

LUMINEX CORPORATION

2021 Annual Meeting
Vote by May 19, 2021
11:59 PM ET



D42140-P53339

You invested in LUMINEX CORPORATION and it's time to vote!

You have the right to vote on proposals being presented at the Annual Meeting. **This is an important notice regarding the availability of proxy material for the stockholder meeting to be held on May 20, 2021.**

Get informed before you vote

View the 2021 Notice and Proxy Statement and Annual Report on Form 10-K online OR you can receive a free paper or email copy of the material(s) by requesting prior to May 6, 2021. If you would like to request a copy of the material(s) for this and/or future stockholder meetings, you may (1) visit www.ProxyVote.com, (2) call 1-800-579-1639 or (3) send an email to sendmaterial@proxyvote.com. If sending an email, please include your control number (indicated below) in the subject line. Unless requested, you will not otherwise receive a paper or email copy.



For complete information and to vote, visit www.ProxyVote.com

Control #

Smartphone users

Point your camera here and
vote without entering a
control number



Vote in Person at the Meeting*

May 20, 2021
10:00 AM CDT

Luminex Corporation
12212 Technology Boulevard
Austin, Texas 78727
See the last page of the Proxy Statement
for directions to the meeting

*Please check the meeting materials for any special requirements for meeting attendance. At the meeting, you will need to request a ballot to vote these shares.

THIS IS NOT A VOTABLE BALLOT

This is an overview of the proposals being presented at the upcoming stockholder meeting. Please follow the instructions on the reverse side to vote these important matters.

| Voting Items | Board Recommends |
|---|---------------------|
| 1. Election of Directors Nominees: | |
| 1a. Edward A. Ogunro, Ph.D. | ☑ For |
| 1b. Kevin M. McNamara | ☑ For |
| 2. Advisory vote to approve named executive officer compensation. | ☑ For |
| 3. Approval of the amendment and restatement of the Luminex Corporation Employee Stock Purchase Plan. | ☑ For |
| 4. Approval of the amendment and restatement of the Luminex Corporation 2018 Equity Incentive Plan. | ☑ For |
| 5. Ratification of the appointment of Ernst & Young LLP as the Company's independent registered public accounting firm for fiscal 2021. | ☑ For |

NOTE: The proxies are further authorized to vote in their discretion on such other business as may properly come before the meeting or any adjournment thereof.

Prefer to receive an email instead? While voting on www.ProxyVote.com, be sure to click "Sign up for E-delivery".