

| OMB APPROVAL | |
|--|-----------|
| OMB Number: | 3235-0287 |
| Estimated average burden hours per response: | 0.5 |

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

| | | | | | | | | | |
|--|---------------|-----------|---|--|--|---|--|--|--|
| 1. Name and Address of Reporting Person* <u>Bennett Todd C.</u> | | | 2. Issuer Name and Ticker or Trading Symbol <u>LUMINEX CORP [LMNX]</u> | | | 5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director _____ 10% Owner _____ Officer (give title below) _____ Other (specify below) _____ <u>SVP, Global Sales & Cust Ops</u> | | | |
| (Last) | (First) | (Middle) | 3. Date of Earliest Transaction (Month/Day/Year) <u>02/25/2021</u> | | | 6. Individual or Joint/Group Filing (Check Applicable Line) <input checked="" type="checkbox"/> Form filed by One Reporting Person Form filed by More than One Reporting Person _____ | | | |
| 12212 TECHNOLOGY BLVD | | | 4. If Amendment, Date of Original Filed (Month/Day/Year) | | | | | | |
| (Street) | <u>AUSTIN</u> | <u>TX</u> | <u>78727</u> | | | | | | |
| (City) | (State) | (Zip) | | | | | | | |

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

| 1. Title of Security (Instr. 3) | 2. Transaction Date (Month/Day/Year) | 2A. Deemed Execution Date, if any (Month/Day/Year) | 3. Transaction Code (Instr. 8) | | 4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5) | | | 5. Amount of Securities Beneficially Owned Following Reported Transaction(s) (Instr. 3 and 4) | 6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) | 7. Nature of Indirect Beneficial Ownership (Instr. 4) |
|---------------------------------|--------------------------------------|--|--------------------------------|---|---|------------|-----------|---|--|---|
| | | | Code | V | Amount | (A) or (D) | Price | | | |
| Common Stock | 02/25/2021 | | M | | 745 | A | \$15.93 | 34,878 | D | |
| Common Stock | 02/25/2021 | | M | | 8,507 | A | \$16.74 | 43,385 | D | |
| Common Stock | 02/25/2021 | | M | | 31,285 | A | \$19.13 | 74,670 | D | |
| Common Stock | 02/25/2021 | | M | | 18,771 | A | \$19.13 | 93,441 | D | |
| Common Stock | 02/25/2021 | | M | | 43,955 | A | \$18.04 | 137,396 | D | |
| Common Stock | 02/25/2021 | | M | | 12,848 | A | \$21.98 | 150,244 | D | |
| Common Stock | 02/25/2021 | | M | | 5,396 | A | \$24.43 | 155,640 | D | |
| Common Stock | 02/25/2021 | | S | | 40,537 ⁽¹⁾ | D | \$33.0702 | 115,103 | D | |
| Common Stock | 02/25/2021 | | S | | 80,970 ⁽¹⁾ | D | \$34.5501 | 34,197 ⁽²⁾ | D | |

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

| 1. Title of Derivative Security (Instr. 3) | 2. Conversion or Exercise Price of Derivative Security | 3. Transaction Date (Month/Day/Year) | 3A. Deemed Execution Date, if any (Month/Day/Year) | 4. Transaction Code (Instr. 8) | | 5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5) | | 6. Date Exercisable and Expiration Date (Month/Day/Year) | | 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4) | | 8. Price of Derivative Security (Instr. 5) | 9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4) | 10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) | 11. Nature of Indirect Beneficial Ownership (Instr. 4) |
|--|--|--------------------------------------|--|--------------------------------|---|--|--------|--|-----------------|---|----------------------------|--|--|---|--|
| | | | | Code | V | (A) | (D) | Date Exercisable | Expiration Date | Title | Amount or Number of Shares | | | | |
| Stock Option (Right To Buy) | \$15.93 | 02/25/2021 | | M | | | 745 | 03/11/2017 | 03/11/2022 | Common Stock | 745 | \$15.93 | 0 | D | |
| Stock Option (Right To Buy) | \$16.74 | 02/25/2021 | | M | | | 8,507 | 07/16/2017 | 10/01/2022 | Common Stock | 8,507 | \$16.74 | 0 | D | |
| Stock Option (Right To Buy) | \$19.13 | 02/25/2021 | | M | | | 31,285 | 03/22/2017 | 03/22/2023 | Common Stock | 31,285 | \$19.13 | 0 | D | |
| Stock Option (Right To Buy) | \$19.13 | 02/25/2021 | | M | | | 18,771 | 03/22/2017 | 03/22/2023 | Common Stock | 18,771 | \$19.13 | 0 | D | |
| Stock Option (Right To Buy) | \$18.04 | 02/25/2021 | | M | | | 43,955 | 03/10/2018 | 03/10/2024 | Common Stock | 43,955 | \$18.04 | 14,652 | D | |
| Stock Option (Right To Buy) | \$21.98 | 02/25/2021 | | M | | | 12,848 | 03/12/2019 | 03/12/2025 | Common Stock | 12,848 | \$21.98 | 12,849 | D | |
| Stock Option (Right To Buy) | \$24.43 | 02/25/2021 | | M | | | 5,396 | 03/12/2020 | 03/12/2026 | Common Stock | 5,396 | \$24.43 | 16,191 | D | |

Explanation of Responses:

1. This sale is being made pursuant to a trading plan intended to comply with Rule 10b5-1 adopted by the seller on August 26, 2020. The representation regarding the seller's knowledge of material adverse non-public information speaks as of the date that the aforementioned trading plan was adopted.

2. This balance includes a total of 64 shares obtained under the Luminex Corporation Employee Stock Purchase Plan. 34 shares obtained March 29, 2020 and 30 shares obtained November 30, 2020.

Remarks:

/S/ HARRISS T.CURRIE AS
ATTORNEY-IN-FACT FOR 02/26/2021
TODD C. BENNETT

** Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

* If the form is filed by more than one reporting person, see Instruction 4 (b)(v).

** Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).

Note: File three copies of this Form, one of which must be manually signed. If space is insufficient, see Instruction 6 for procedure.

Persons who respond to the collection of information contained in this form are not required to respond unless the form displays a currently valid OMB Number.

POWER OF ATTORNEY

Know all by these presents, that the undersigned hereby constitutes and appoints Harriss T. Currie and David S. Reiter, signing singly, the undersigned's true and lawful attorney-in-fact to:

- (1) execute for and on behalf of the undersigned, in the undersigned's capacity as an officer and/or director of Luminex Corporation (the "Company"), Forms 3, 4, and 5 in accordance with Section 16(a) of the Securities Exchange Act of 1934 and the rules thereunder;
- (2) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such Form 3, 4, or 5, complete and execute any amendment or amendments thereto, and file such form with the United States Securities and Exchange Commission and any stock exchange or similar authority; and
- (3) take any other action of any type whatsoever in connection with the foregoing which, in the opinion of such attorney-in-fact, may be of benefit to, in the best interest of, or legally required of, the undersigned, it being understood that the documents executed by such attorney-in-fact on behalf of the undersigned pursuant to this Power of Attorney shall be in such form and shall contain such terms and conditions as such attorney-in-fact may approve in such attorney-in-fact's discretion.

The undersigned hereby grants to each such attorney-in-fact full power and authority to do and perform any and every act and thing whatsoever requisite, necessary, or proper to be done in the exercise of any of the rights and powers herein granted, as fully to all intents and purposes as the undersigned might or could do if personally present, with full power of substitution or revocation, hereby ratifying and confirming all that such attorney-in-fact, or such attorney-in-fact's substitute or substitutes, shall lawfully do or cause to be done by virtue of this power of attorney and the rights and powers herein granted. The undersigned acknowledges that the foregoing attorneys-in-fact, in serving in such capacity at the request of the undersigned, are not assuming, nor is the Company assuming, any of the undersigned's responsibilities to comply with Section 16 of the Securities Exchange Act of 1934.

This Power of Attorney shall remain in full force and effect until the undersigned is no longer required to file Forms 3, 4, and 5 with respect to the undersigned's holdings of and transactions in securities issued by the Company, unless earlier revoked by the undersigned in a signed writing delivered to the foregoing attorneys-in-fact.

IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed as of this 15th day of November, 2016.

/S/ Todd C. Bennett

Name: Todd C. Bennett