

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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

| | | |
|---|---|---|
| 1. Name and Address of Reporting Person* <u>LOEWENBAUM G WALTER II</u> (Last) (First) (Middle) 12212 TECHNOLOGY BOULEVARD (Street) AUSTIN TX 78727 (City) (State) (Zip) | 2. Issuer Name and Ticker or Trading Symbol <u>LUMINEX CORP [LMNX]</u> | 5. Relationship of Reporting Person(s) to Issuer (Check all applicable) <input checked="" type="checkbox"/> Director 10% Owner Officer (give title below) Other (specify below) |
| | 3. Date of Earliest Transaction (Month/Day/Year) 07/01/2020 | |
| 4. If Amendment, Date of Original Filed (Month/Day/Year) | | |

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 | | | | | | | | 550,189 | D | |
| Common Stock | 07/01/2020 | 07/01/2020 | s | | 7,853 ⁽¹⁾ | D | \$32.8847 ⁽²⁾ | 801,562 | I | Spouse ⁽³⁾ |
| Common Stock | 07/02/2020 | 07/02/2020 | s | | 15,800 ⁽¹⁾ | D | \$32.8577 ⁽⁴⁾ | 785,762 | I | Spouse ⁽³⁾ |
| Common Stock | | | | | | | | 96,614 | I | Partnership ⁽³⁾ |
| Common Stock | | | | | | | | 428,520 | I | Trusts ⁽³⁾ |

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) |
|--|--|--------------------------------------|--|--------------------------------|--|-----|--|-----------------|---|--|--|---|--|
| | | | | | (A) | (D) | Date Exercisable | Expiration Date | | | | | |

Explanation of Responses:

- This sale is being made pursuant to a trading plan intended to comply with Rule 10b5-1, adopted by the seller on May 27, 2020 for shares acquired as founder's stock prior to the IPO on March 30, 2000. 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.
- The price reported in Column 4 is a weighted average price. These shares were sold in multiple transactions at prices ranging from \$32.85 to \$32.925 inclusive. Mr. Loewenbaum undertakes to provide to Luminex Corporation, any security holder of Luminex Corporation, or the staff of the Securities and Exchange Commission, upon request, full information regarding the number of shares sold at each separate price within the range set forth in this footnote to this Form 4.
- The reporting person disclaims beneficial ownership of these securities except to the extent of his pecuniary interest therein, and the inclusion of these shares in this report shall not be deemed an admission of beneficial ownership of all of the reported shares for purposes of Section 16 or for any other purpose.
- The price reported in Column 4 is a weighted average price. These shares were sold in multiple transactions at prices ranging from \$32.85 to \$32.90 inclusive. Mr. Loewenbaum undertakes to provide to Luminex Corporation, any security holder of Luminex Corporation, or the staff of the Securities and Exchange Commission, upon request, full information regarding the number of shares sold at each separate price within the range set forth in this footnote to this Form 4.

Remarks:

/S/ HARRISS T. CURRIE, AS
ATTORNEY-IN-FACT FOR 07/02/2020
G. WALTER
LOEWENBAUM II

** 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.