

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>Rew Richard W. II</u> (Last) (First) (Middle) <u>12212 TECHNOLOGY BLVD</u> (Street) <u>AUSTIN TX 78727</u> (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) Director 10% Owner <input checked="" type="checkbox"/> Officer (give title below) Other (specify below) <u>SVP, Gen Counsel & Corp Secy</u>
	3. Date of Earliest Transaction (Month/Day/Year) <u>03/11/2021</u>	
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 Beneficial Ownership (Instr. 4)
			Code	V	Amount	(A) or (D)	Price			
Common Stock	03/11/2021		F		295 ⁽¹⁾	D	\$32.45	32,858	D	
Common Stock	03/12/2021		F		965 ⁽²⁾	D	\$32.76	31,893	D	
Common Stock	03/12/2021		A		14,957	A	\$32.76	47,272 ⁽³⁾	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					
Stock Option (Right To Buy)	\$23.09	03/12/2021		A		66,490 ⁽⁴⁾		03/12/2021	03/12/2027	Common Stock	66,490	66,490	D	

Explanation of Responses:

- This disposition relates to withholding for taxes on 1,213 vesting shares associated with a previously reported issuance.
- This disposition relates to withholding for taxes on 6,188 vesting shares associated with a previously reported issuance.
- This balance includes a total of 298 shares obtained November 30, 2020 under the Luminex Corporation Employee Stock Purchase Plan, and 124 shares obtained under a personal dividend reinvestment plan.
- The Board of Directors of Luminex Corporation determined the level of performance and number of shares earned for this performance option. The option will vest annually in four equal installments beginning 03/12/21.

Remarks:

/S/ HARRISS T.CURRIE AS
ATTORNEY-IN-FACT FOR 03/15/2021
RICHARD W. REW 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.