

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>Rew Richard W. II</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) <input checked="" type="checkbox"/> Director <input type="checkbox"/> Officer (give title below) <input type="checkbox"/> 10% Owner <input type="checkbox"/> Other (specify below) <u>SVP, Gen Counsel & Corp Secy</u>
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year) <u>07/17/2020</u>		
12212 TECHNOLOGY BLVD			4. If Amendment, Date of Original Filed (Month/Day/Year)		
(Street)	<u>AUSTIN</u>	<u>TX</u>	<u>78727</u>	6. Individual or Joint/Group Filing (Check Applicable Line) <input checked="" type="checkbox"/> Form filed by One Reporting Person <input type="checkbox"/> Form filed by More than One Reporting Person	
(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	07/17/2020		M		11,175	A	\$15.93	43,984	D	
Common Stock	07/17/2020		S		11,175 ⁽¹⁾	D	\$36.93	33,153 ⁽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						Amount or Number of Shares
Stock Option (Right To Buy)	\$15.93	07/17/2020		M			11,175	03/11/2017	03/11/2022	Common Stock	11,175	\$15.93	33,527	D	

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 22, 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.
- This balance includes a total of 344 shares obtained under the Luminex Corporation Employee Stock Purchase Plan on May 29, 2020.

Remarks:

/S/ HARRISS T.CURRIE AS
ATTORNEY-IN-FACT FOR 07/17/2020
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.