

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>Fairchild Nancy</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) _____ SVP, Human Resources		
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year) <u>07/06/2020</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		
<u>12212 TECHNOLOGY BLVD</u>			4. If Amendment, Date of Original Filed (Month/Day/Year)					
(Street)	<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	07/06/2020		M		3,475	A	\$19.25	60,899	D	
Common Stock	07/06/2020		S		3,475 ⁽¹⁾	D	\$33	57,424	D	
Common Stock	07/06/2020		M		20,000	A	\$15.93	77,424	D	
Common Stock	07/06/2020		S		20,000 ⁽¹⁾	D	\$33	58,111 ⁽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
Stock Option (Right To Buy)	\$19.25	07/06/2020		M			3,475	08/14/2013	10/01/2022	Common Stock	3,475	\$0	0	D	
Stock Option (Right To Buy)	\$15.93	07/06/2020		M			20,000	03/11/2017	03/11/2022	Common Stock	20,000	\$0	32,153	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 20, 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 687 shares obtained under the Luminex Corporation Employee Stock Purchase Plan on May 29, 2020.

Remarks:

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

** 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.