

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 _____ <input checked="" type="checkbox"/> Officer (give title below) _____ Other (specify below) _____ <u>SVP, Human Resources</u>	
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year) <u>07/15/2020</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 _____	
12212 TECHNOLOGY BLVD			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	07/15/2020		M		18,399	A	\$19.13	76,510	D	
Common Stock	07/15/2020		S		18,399 ⁽¹⁾	D	\$36.0087	58,111	D	
Common Stock	07/15/2020		M		32,153	A	\$15.93	90,264	D	
Common Stock	07/15/2020		S		32,153 ⁽¹⁾	D	\$36.0056	58,111	D	
Common Stock	07/15/2020		M		5,973	A	\$17.24	64,084	D	
Common Stock	07/15/2020		S		5,973 ⁽¹⁾	D	\$36	58,111	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)	\$19.13	07/15/2020		M			18,399	03/22/2017	03/22/2023	Common Stock 18,399	\$19.13	372	D	
Stock Option (Right To Buy)	\$15.93	07/15/2020		M			32,153	03/11/2017	03/11/2022	Common Stock 32,153	\$15.93	0	D	
Stock Option (Right To Buy)	\$17.24	07/15/2020		M			5,973	03/08/2014	03/08/2023	Common Stock 5,973	\$0	0	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 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.

Remarks:

/S/ HARRISS T.CURRIE AS
ATTORNEY-IN-FACT FOR 07/16/2020
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.