

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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>Myers Randall</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, Global Mfg and Quality</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 Indirect Beneficial Ownership (Instr. 4)
			Code	V	Amount	(A) or (D)	Price			
Common Stock	03/11/2021		F		295 ⁽¹⁾	D	\$32.45	33,843	D	
Common Stock	03/12/2021		F		1,596 ⁽²⁾	D	\$32.76	32,247	D	
Common Stock	03/12/2021		A		16,788	A	\$32.76	49,511 ⁽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	Title	Amount or Number of Shares				
Stock Option (Right To Buy)	\$23.09	03/12/2021		A		72,976 ⁽⁴⁾		03/12/2021	03/12/2027	Common Stock	72,976	\$23.09	72,976	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 6,558 vesting shares associated with a previously reported issuance.
- This balance includes 476 shares obtained November 30, 2020 under the Luminex Corporation Employee Stock Purchase 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/ Harris T. Currie As Attorney-in-Fact 03/15/2021
for Randall Myers

** 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.

POWER OF ATTORNEY

Know all by these presents, that the undersigned hereby constitutes and appoints Harriss T. Currie and Richard W. Rew II, signing singly
(1) execute for and on behalf of the undersigned, in the undersigned's capacity as an officer and/or director of Luminex Corporation (the "Company")
(2) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such
(3) take any other action of any type whatsoever in connection with the foregoing which, in the opinion of such attorney-in-fact, may be of benefit to the undersigned.

The undersigned hereby grants to each such attorney-in-fact full power and authority to do and perform any and every act and thing whatsoever that the undersigned may lawfully do or perform.

This Power of Attorney shall remain in full force and effect until the undersigned is no longer required to file Forms 3, 4, and 5 with the SEC.
IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed as of March 16, 2015.

/S/ RANDALL MYERS

Name: Randall Myers