

**UNITED STATES
SECURITIES AND EXCHANGE COMMISSION**

Washington, D.C. 20549

FORM 8-K

**CURRENT REPORT
Pursuant to Section 13 OR 15(d)
of The Securities Exchange Act of 1934**

Date of Report (Date of earliest event reported): May 24, 2021

Luminex

LUMINEX CORPORATION

(Exact name of registrant as specified in its charter)

Delaware
(State or other jurisdiction
of incorporation)

000-30109
(Commission
File Number)

74-2747608
(IRS Employer
Identification No.)

12212 Technology Blvd., Austin, Texas
(Address of principal executive offices)

78727
(Zip Code)

(512) 219-8020

Registrant's Telephone Number, Including Area Code

None

(Former name or former address, if changed since last report.)

Check the appropriate box below if the Form 8-K filing is intended to simultaneously satisfy the filing obligation of the registrant under any of the following provisions:

- Written communications pursuant to Rule 425 under the Securities Act (17 CFR 230.425)
- Soliciting material pursuant to Rule 14a-12 under the Exchange Act (17 CFR 240.14a-12)
- Pre-commencement communications pursuant to Rule 14d-2(b) under the Exchange Act (17 CFR 240.14d-2(b))
- Pre-commencement communications pursuant to Rule 13e-4(c) under the Exchange Act (17 CFR 240.13e-4(c))

Securities registered pursuant to Section 12(b) of the Act:

Title of each class	Trading Symbol(s)	Name of exchange on which registered
Common Stock, \$0.001 par value	LMNX	The Nasdaq Global Select Market

Indicate by check mark whether the registrant is an emerging growth company as defined in Rule 405 of the Securities Act of 1933 (§230.405 of this chapter) or Rule 12b-2 of the Exchange Act (§240.12b-2 of this chapter).

Emerging growth company

If an emerging growth company, indicate by check mark if the registrant has elected not to use the extended transition period for complying with any new or revised financial accounting standards provided pursuant to Section 13(a) of the Exchange Act.

Item 8.01. Other Events.

On May 24, 2021, the waiting period under the Hart-Scott-Rodino Antitrust Improvements Act of 1976, as amended (the “HSR Act”), expired in connection with the previously announced proposed merger (the “Merger”) of Diagonal Subsidiary Inc., a Delaware corporation and wholly owned subsidiary of DiaSorin S.p.A., a *società per azioni* organized under the laws of the Republic of Italy (“Parent”), with and into Luminex Corporation, a Delaware corporation (the “Company”), with the Company surviving the Merger as a wholly owned subsidiary of Parent.

The expiration of the waiting period under the HSR Act satisfies one of the conditions to the closing of the Merger. The closing of the Merger remains subject to the satisfaction or waiver of the remaining conditions to the Merger set forth in the related Agreement and Plan of Merger (the “Merger Agreement”), dated as of April 11, 2021, by and among the Company, Parent and Merger Sub, including adoption of the Merger Agreement by the Company’s stockholders and the receipt of certain other regulatory approvals.

Forward-looking statements

This Current Report on Form 8-K, and the documents to which the Company refers you in this communication, contains not only historical information, but also forward-looking statements made pursuant to the safe-harbor provisions of the Private Securities Litigation Reform Act of 1995. These forward-looking statements represent the Company’s expectations or beliefs concerning future events, including the timing of the transaction and other information relating to the transaction. Forward-looking statements include information concerning possible or assumed future results of operations of the Company, the expected completion and timing of the transaction and other information relating to the transaction. Without limiting the foregoing, the words “believes,” “anticipates,” “plans,” “expects,” “intends,” “forecasts,” “should,” “estimates,” “contemplate,” “future,” “goal,” “potential,” “predict,” “project,” “projection,” “may,” “will,” “could,” “should,” “would,” “assuming” and similar expressions are intended to identify forward-looking statements. You should read statements that contain these words carefully. They discuss the Company’s future expectations or state other forward-looking information and may involve known and unknown risks over which the Company has no control. Those risks include, (i) the risk that the transaction may not be completed in a timely manner or at all, which may adversely affect the Company’s business and the price of the common stock of the Company, (ii) the failure to satisfy the conditions to the consummation of the transaction, including the adoption of the Merger Agreement by the stockholders of the Company and the receipt of regulatory approvals from various governmental entities (including any conditions, limitations or restrictions placed on these approvals) and the risk that one or more governmental entities may deny approval, (iii) the occurrence of any event, change or other circumstance that could give rise to the termination of the Merger Agreement, (iv) the risk that the definitive Merger Agreement may be terminated in circumstances that require the Company to pay a termination fee; (v) risks regarding the failure to obtain the necessary financing to complete the Merger, (vi) the effect of the announcement or pendency of the transaction on the Company’s business relationships, operating results and business generally, (vii) risks that the proposed transaction disrupts current plans and operations, (viii) risks related to diverting management’s attention from the Company’s ongoing business operations, and (ix) the outcome of any legal proceedings that may be instituted against the Company related to the Merger Agreement or the transaction. Forward-looking statements speak only as of the date of this communication or the date of any document incorporated by reference in this document. Further risks that could cause actual results to differ materially from those matters expressed in or implied by such forward-looking statements are described in the Company’s Securities and Exchange Commission (the “SEC”) reports, including but not limited to the risks described in the Company’s Annual Report on Form 10-K for its fiscal year ended December 31, 2020. Except as required by applicable law or regulation, the Company does not undertake to update these forward-looking statements to reflect future events or circumstances.

Additional Information and Where to Find It

On May 20, 2021, in connection with the Merger, the Company prepared and filed with the SEC a definitive proxy statement for the Company’s stockholders in connection with the transactions contemplated by the Merger Agreement. The proxy statement has also been mailed to the Company’s stockholders of record on the record date. INVESTORS AND SECURITY HOLDERS OF THE COMPANY ARE URGED TO READ THESE MATERIALS (INCLUDING ANY AMENDMENTS OR SUPPLEMENTS THERETO) AND ANY OTHER RELEVANT DOCUMENTS IN CONNECTION WITH THE MERGER THAT THE COMPANY WILL FILE WITH THE SEC WHEN THEY BECOME AVAILABLE BECAUSE THEY WILL CONTAIN IMPORTANT INFORMATION ABOUT THE COMPANY AND THE MERGER. The definitive proxy statement, the preliminary proxy statement and other relevant materials in connection with the Merger, and any other

documents filed by the Company with the SEC, may be obtained free of charge at the SEC's website (<http://www.sec.gov>) or at the Company's website luminexcorp.com or by writing to the Company's Secretary at 12212 Technology Blvd., Austin, Texas 78727.

Participants in the Solicitation

The Company and its directors and executive officers may be deemed to be participants in the solicitation of proxies from the Company's stockholders with respect to the Merger. Information about the Company's directors and executive officers and their ownership of the Company's common stock is set forth in the proxy statement on Schedule 14A filed with the SEC on March 31, 2021 and the Company's Annual Report on Form 10-K for the fiscal year ended December 31, 2020. To the extent that such individual's holdings of the Company's common stock have changed since the amounts printed in the Company's proxy statement, such changes have been or will be reflected on Statements of Change in Ownership on Form 4 filed with the SEC. Information regarding the identity of the potential participants, and their direct or indirect interests in the Merger, by security holdings or otherwise, are set forth in the proxy statement and other materials filed with SEC in connection with the Merger.

SIGNATURES

Pursuant to the requirements of the Securities Exchange Act of 1934, as amended, the registrant has duly caused this report to be signed on its behalf by the undersigned hereunto duly authorized.

LUMINEX CORPORATION

By: /s/ Harriss T. Currie

Harriss T. Currie

Chief Financial Officer, Senior Vice President of
Finance (Principal Financial Officer)

Date: May 26, 2021