

The information contained herein is being provided pursuant to the requirements of Section 6045B of the Internal Revenue Code of 1986, as amended. The information in this document does not constitute tax advice and should not be construed to take into account any shareholder's specific circumstances. Holders and nominees should consult their own tax advisors regarding the particular tax consequences of the organizational action (as described in this document) to them, including the applicability and effect of all U.S. federal, state, and local and foreign tax laws.

Luminex Corporation
Attachment to Internal Revenue Service Form 8937

On May 7, 2020, Luminex Corporation (the “**Company**”) issued 3.00% convertible senior notes due 2025 (the “**Notes**”). The Company has provided a notice of adjustment to the conversion rate applicable to the Notes.

Section 6045B of the Internal Revenue Code and the related Treasury Regulations provide that effective January 1, 2016, if a company pays a cash dividend with respect to stock resulting in a conversion rate adjustment on a convertible debt instrument that is treated as a deemed distribution under 305(b)(2) and (c) of the Code, an information return on Form 8937 must be filed with the Internal Revenue Service or, in lieu of filing, posted to the company’s public website.

The information in this document does not constitute tax advice and is not intended or written to be used, and cannot be used, for the purpose of avoiding penalties under the Code. Holders of the Notes should consult their own tax advisors regarding the particular tax consequences of the conversion rate adjustment to them, including the applicability and effect of all U.S. federal, state and local and foreign tax laws. Such information is not intended to be a complete analysis or description of all potential federal or other tax consequences of the conversion rate adjustment.

Form 8937, Part II, Line 14

Effective immediately after the open of business on December 22, 2020 (i.e., the ex-dividend date of the dividend paid on January 14, 2021), in connection with the cash dividends previously announced by the Company on December 3, 2020, the conversion rate of the Notes was adjusted from 22.8918 to 22.9009 shares of Company’s common stock per \$1,000 principal amount of the Notes.

Form 8937, Part II, Line 15

In connection with the Company’s payment of the cash dividend described in Line 14, the tax basis adjustment is an increase of \$0.2310 per \$1,000 of principal amount of Notes. The Company expects that it will have sufficient current or accumulated earnings and profits. Accordingly, as a result of the increase in the conversion rate of the Notes, each holder of the Notes is expected to include \$0.2310 in gross income as a dividend per \$1,000 principal amount of the Notes with a resulting basis increase of the same amount.

Form 8937, Part II, Line 16

The conversion rate adjustment is calculated based on the following formula where such adjustment would require certain conditions to be met to make it effective as further specified in article 14.04(d) of the Notes' indenture:

$$CR_1 = CR_0 \times \frac{SP_0 - T}{SP_0 - C}$$

$$\begin{aligned} \text{The basis adjustment} &= SP_0 * (CR_1 - CR_0) \\ &= \$25.3900 * (22.9009 - 22.8918) \\ &= \$0.2310 \text{ per } \$1,000 \text{ principal amount of the Notes} \end{aligned}$$

Accordingly, for illustrative purposes, a holder of the Notes with a principal amount of \$1,000 is expected to increase its basis in the Notes by \$0.2310 as a result of the inclusion of the dividend in gross income for U.S. federal income tax purposes.

Form 8937, Part II, Line 17

Sections 305(b)(2), 305(c), 301(c) and 301(d).

Form 8937, Part II, Line 18

There is no associated loss recognition.

Form 8937, Part II, Line 19

The reportable tax year for this conversion rate adjustment is 2020 for a calendar year taxpayer.