

**CHARTER OF THE
EXECUTIVE COMMITTEE
OF THE BOARD OF DIRECTORS
OF
LUMINEX CORPORATION**

Purpose

The purpose of the Executive Committee (the “Committee”) of the Board of Directors (the “Board”) of Luminex Corporation (the “Company”) shall be to exercise the power and authority of the Board between meetings or otherwise when it is not reasonably possible to convene a special meeting of the full Board or when the Board has delegated a specific matter to it for action, subject to the limitations provided herein. Notwithstanding anything herein to the contrary, the Board may limit or qualify the powers of the Committee at any time.

Powers and Duties

1. The Committee shall exercise all of the powers and authority of the full Board between meetings or otherwise when it is not reasonably possible to convene a special meeting of the full Board or when the Board has delegated a specific matter to it for action, subject to such limitations as the Board and/or applicable law may from time to time impose; provided that in no event does the Committee have the power or authority to (i) declare dividends or other corporate distributions; (ii) remove members of the Board (iii) amend the Company’s Certificate of Incorporation or Bylaws; (iv) adopt an agreement or plan of merger, share exchange, or consolidation to which the Company is a party; or (v) recommend to the stockholders of the Company any action which requires stockholder approval.
2. Review and evaluate certain transactions or special projects as requested by the Board and, if so requested, bring recommendations to the Board regarding those transactions or projects.
3. The Committee shall have the authority to obtain advice and assistance from internal and external legal, accounting or other consultants or advisors it may deem necessary or appropriate in connection with carrying out its duties and to approve the fees and other retention terms related to any such external counsel, consultants and advisors.
4. The Committee may form and delegate authority to subcommittees, of not less than two Board members, where appropriate.
5. The Committee shall periodically review and reassess the adequacy of this Charter and recommend any proposed changes to the Board for approval. At least annually the Committee shall evaluate its own performance.

Composition of Committee

The Committee shall be appointed by the Board and shall consist of not less than three directors, including a Committee Chairman. Directors will serve on the Committee at the pleasure of the Board and for such terms as the Board may determine.

Meetings

The Committee shall meet as frequently as necessary or appropriate. Special meetings of the Committee may be called on two hours notice by any member of the Committee. A majority of the Committee shall constitute a quorum, and the Committee shall act only on the affirmative vote of a majority of the members present at the meeting. The Committee shall keep such records of its activities as it shall deem appropriate. A summary of the actions taken by the Committee shall be presented to the Board in the regular course of the Board's activities.

Committee Access and Resources

In carrying out its responsibilities, the Committee shall have access to all the Company's books, records, directors, officers, employees, consultants and other advisors.

Dated September 9, 2009