

**Research Travel
Grants****PLEASE READ THESE GUIDELINES
COMPLETELY BEFORE COMPLETING AND
SUBMITTING AN APPLICATION FORM.**

The Luminex Research Travel Grant (RTG) program is intended to defray travel and related expenses incurred by Luminex customers related to customer presentations of xMAP®-related research. Possible presentation opportunities include tradeshow, regional meetings, or Luminex- or partner- hosted seminars.

Benefits to the xMAP researcher are 1) direct communication with Luminex and/or partners who provide hardware and reagents, and 2) funding for travel to national/international society meetings or regional user groups.

Benefits to Luminex and/or partners are 1) direct communication with researchers using xMAP technology, 2) technology credibility through 3rd-party research and evaluations, and 3) exposure of technology and corporate brand to targeted markets.

Definitions.

“Sponsoring Institution” means the employer or sponsor of a researcher where such employer or sponsor is a customer of Luminex.

“Applicant” means an individual who is or has conducted research related to Luminex xMAP Technologies and is designated on a completed RTG Application as a presenter or researcher in a topic of interest to Luminex.

“RTG Application Form” means the form of application attached to these guidelines.

“RTG Program” means the Luminex Research Travel Grant Program set forth in these guidelines.

**General RTG
Program
Guidelines**

Luminex will support key researchers in presenting approved xMAP-related research data, and will work with partners to support such presentations. Guidelines of these collaborations are as follows.

- Travel Grants are subject to review and approval by Luminex in its sole discretion. Applicants must submit a complete and fully executed RTG Application Form, including a proposed abstract, prior to committing to an event. The proposed travel and event must be consistent with these guidelines. Details and instructions for completing and submitting the RTG Application Form are set forth below.
- Luminex will review the application and will promptly advise the Applicant whether the application is approved, declined or whether more information is requested. See details below.
- Once an application is approved, upon completion of an approved presentation, Luminex will reimburse the Applicant for approved expenses.
- Except as otherwise specified in the approval referenced above, Luminex will reimburse up to \$1500 of travel expenses resulting from approved presentations, including the following types of expenses: registration costs, airfare, hotel, meals, or other incidentals during the trip. At the request of the Applicant, Luminex will also work with our partners to cover additional costs associated with approved presentations.

**Acceptable
Research
Topics**

Applications that identify research in the following areas are most likely to be approved for Grants under the RTG Program:

- Novel applications of xMAP technology within the Life Sciences, including Clinical Research and Diagnostics, Drug Discovery and Development, Agriculture, and BioDefense
- Clinically important findings regarding disease mechanisms, diagnosis, or treatment
- Comparative workflow analyses between xMAP-based systems and traditional methods such as ELISA, Western Blot, genetic sequencing, real-time PCR, or microarray
- Correlation studies between accepted platforms such as ELISA or Mass Spectrometry and xMAP technology
- Overview and benefits of xMAP multiplexing versus traditional single-plex platforms or competitive, novel multiplex platforms.

**Travel Grant
Application and
Approval
Process**

Luminex will accept applications for Research Travel Grants (RTG's) from customer researchers and will work with the partner to fund such applications. Here is the approval process:

1. To apply for a Luminex Research Travel Grant, an Applicant must submit a proposed abstract to the Luminex RTG Program Manager using the attached RTG application form. The completed RTG form must include at a minimum the name of the event, the general audience of the event, the general importance of research findings, the duration of the event and the expected travel, lodging and related expenses. The application form must be signed by the Applicant and by an authorized representative of the Sponsoring Institution before submission to Luminex.
2. When Luminex receives a RTG Application Form, the Luminex RTG program manager will evaluate the proposal and mark it as "Approved," "Declined" or "More Information Requested" and return it to the Applicant. If Luminex determines to award a travel grant under this Program based on a completed RTG Application Form, Luminex will sign the final, approved RTG Application Form indicating the maximum amount to be reimbursed by Luminex for approved expenses and return it to the applicant.
3. After the event occurs, the applicant must send Luminex a copy of any written or printed materials used in connection with the presentation together with copies of the direct expenses (invoices, etc.), and proof of payment (copies of checks or charge receipts). Luminex will reimburse applicants for approved expenses up to the amount authorized in the approved RTG Application Form within 30 days following the end of the calendar quarter in which Luminex approves such expenses.

**RTG Program
Conditions**

The Luminex Research Travel Grant Program has the following conditions:

1. This program is for domestic US and Canadian travel only. Separate programs may apply for International partners and distributors.
2. Luminex must pre-approve all travel grant proposals. Luminex reserves the right to not reimburse any expenses incurred before such approval is granted in writing. The researcher has the risk that expenses incurred until formal approval may not be reimbursed.
3. In order to qualify for a Travel Grant, an xMAP Researcher's Sponsoring Institution must be current in its accounts with Luminex to receive any Research Travel Grant funds.
4. A Sponsoring Institution may not directly allocate travel grant funds to outstanding balances with Luminex.
5. Payment is limited to the funds available for the approved travel event and dates. It is not possible to carry over fund requests from one event into the next if funds are leftover from a previous event.
6. Luminex retains all rights in interpreting the rules for this program. Luminex is the final authority, with sole discretion, in making decisions on the implementation of this program and the approval of projects.
7. Luminex may cancel this program at any time with no notice.
8. Submission of projects for reimbursement constitutes agreement to the program and these conditions.

Research Travel Grant Application Form

Research Travel Grant #: _____ *
**(to be assigned by Luminex)*

Name of tradeshow, meeting, or other event:

General audience:

Abstract (include title and use guidelines set by the intended event; (attach additional pages if necessary):

General importance of research findings:

Duration of event:

Expected travel, lodging, and other costs:

Airfare:

Ground Transportation:

Lodging:

Registration Fees:

Meals

Other (please specify):

Date approval by Luminex required

Approval:

Both parties agree to the stated amount of reimbursement for this Travel Grant, agree to the expiration date below, and agree to the general terms and guidelines stated within this document.

(up to \$1500) _____
(Funded Amount by Luminex)

12/31/06 _____
(Program Expiration Date)

Authorization Signatures

Grant Gibson, Ph.D., Luminex RTG Program Manager

<applicant signature>, <company/institution>