## Luminex

# Let's Partner

Unlock new opportunities for your business with Luminex's cutting-edge xMAP<sup>®</sup> Technology—the world's leading multiplex platform. Whether you're looking to commercialize a multiplex assay or transform your existing immunoassay or nucleic acid assay, we offer a flexible, proven, and widely adopted platform to meet your needs.

Embrace efficiency and cost-effectiveness as you leverage Luminex's proprietary xMAP Technology through our innovative licensing model. By partnering with us, you can accelerate your commercial success in a broad spectrum of life science, biopharma, and clinical applications. Let our platform empower your endeavors and open doors to a world of possibilities.

Get ready to take your business to new heights with Luminex—the gateway to a brighter future in the realm of multiplex assays.

#### **Partner Facts**

- >70 Luminex partners globally
- Our partners' kits and testing services span protein and nucleic acid applications for IVD, applied markets, biopharma and research use
- Our partners include industry leaders such as Thermo Fisher Scientific®, MilliporeSigma™, Bio-Techne™, and Bio-Rad®

### Partner Collaboration

- Team up with LuminexPLORE Lab to achieve your custom assay development goals
- Partner with Luminex to leverage your sales and marketing initiatives
- Participate within a scientific community reinforced by the widely validated benefits of this established platform

#### xMAP Benefits Include:

- Tests up to **500 analytes** simultaneously from one sample, in one well
- Proven technology unrivaled in the level of end-user adoption and validation
- Supports a wide range of research, translational, and diagnostic applications with **four instrument options** and a broad install-base
- Yields more data while expending less in time, sample, and cost
- Flexible, reproducible platform with high-throughput and enhanced statistical power
- Efficient, open architecture design enabling the use of commercial kits and internally developed assays
- >100 FDA 510(k) cleared kits
- **>75,000 scientific publications** cite xMAP Technology and/or our partners' products



#### To learn more, visit info.diasorin.com/en-us/research/become-a-partner

©2023 Luminex Corporation. A DiaSorin Company. All rights reserved. Luminex and xMAP are trademarks of Luminex Corporation, registered in the US and other countries. All other trademarks are the property of their respective owners.