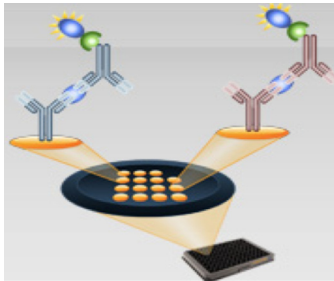


Multiplex ELISA

Q-Plex™ Multiplex Cytokine Arrays

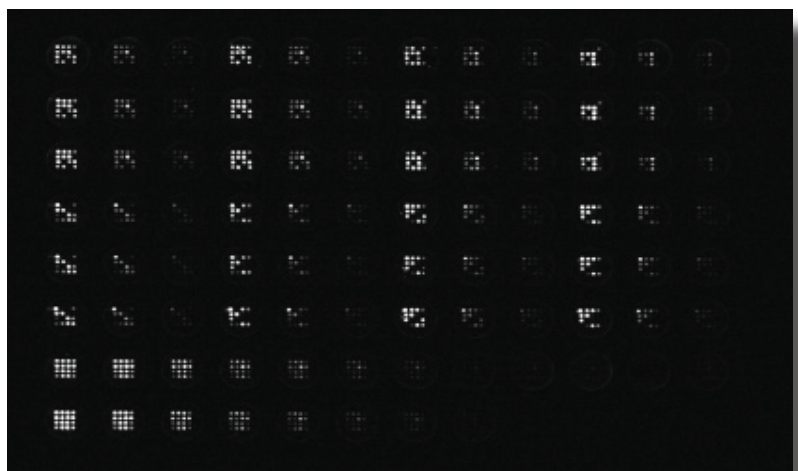
Effective Cytokine Profiling For Less Cost



Product Features:

- Up to 16 cytokines can be assayed in 2.5 hours
- Detection ranges from 30 pg/ml to less than 1 pg/ml
- Less than 30 µl sample requirement
- Compatible with serum, plasma, cell culture supernatant, tissue homogenates, lysates, nasal lavage, tears and urine

PBL InterferonSource is pleased to be the worldwide distributor of Q-Plex™ Multiplex Arrays by Quansys Biosciences. The Q-Plex™ (multiplex ELISA) arrays are fully quantitative ELISA-based tests where up to 16 distinct capture antibodies have been absorbed to each well of a 96-well plate in a defined array. Each spot represents a different distinct capture antibody population. The ability to assess multiple cytokines in one assay means more time to generate cohesive information and knowledge for researchers. Multiplex has consistently proven to be more efficient than traditional ELISAs in terms of time, cost and sample amount. However, not all multiplexing methods are equal. The Q-Plex™ Multiplex Arrays are applied in 96-well plates at a cost of under \$1.00 per data point, whereas other technologies employ membranes, chips, or even bead-based arrays at a cost that is two- to four-times higher per data point. The Quansys Q-Plex™ Array is an invaluable tool for the simultaneous detection of multiple analytes. It provides a platform that overcomes the shortfalls of both conventional ELISA assays and predecessor multiplex technologies to provide the researcher an easy and cost-effective means of exploring the complexity and range of many biological analytes. In short, the Quansys Q-Plex™ Array is faster, simpler and far less expensive than traditional ELISAs as well as other multiplex technologies and yields much more data using only minimal sample amounts. For more information, please visit our web site at: www.interferonsource.com/multiplex or contact us at: info@interferonsource.com.



Example Plate

Q-Plex™ Multiplex Arrays

Catalog No.	Product Description	Antigen
110351HU	Q-Plex™ Human Cytokine-Stripwells™ (16-plex)	IL-1 α , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, IL-15, IL-17, IL-23, IFN- γ , TNF- α , TNF- β
110451HU	Q-Plex™ Human Cytokine-Inflammation (9-plex)	IL-1 α , IL-1 β , IL-2, IL-4, IL-6, IL-8, IL-10, IFN- γ , TNF- α
110951HU	Q-Plex™ Human Cytokine Screen (16-plex)	IL-1 α , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, IL-15, IL-17, IL-23, IFN- γ , TNF- α , TNF- β
130251HU	Q-Plex™ Human Cytokine-IR (9-plex, infrared)	IL-1 α , IL-1 β , IL-2, IL-4, IL-6, IL-10, IL-12p70, IFN- γ , TNF- α
130951HU	Q-Plex™ Human Cytokine-Screen IR (16-plex)	IL-1 α , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, IL-15, IL-17, IFN- γ , MCP-1, TNF- α , TNF- β
150251HU	Q-Plex™ Human Angiogenesis (9-plex)	Ang-2, FGF basic, HGF, IL-8, PDGF-BB, TIMP-1, TIMP-2, TNF- α , VEGF
110951MS	Q-Plex™ Mouse Cytokine-Screen (16-plex)	IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-10, IL-12p70, IL-17, MCP-1, IFN- γ , TNF- α , MIP-1 α , GM-CSF, RANTES
110451MS	Q-Plex™ Mouse Cytokine-Inflammation (14-plex)	IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-6, IL-10, IL-12p70, IL-17, MCP-1, TNF- α , MIP-1 α , GM-CSF, RANTES
130251MS	Q-Plex™ Mouse Cytokine-IR (9-plex, Infrared)	IL-1 α , IL-1 β , IL-2, IL-4, IL-6, IL-10, IL-12p70, IFN- γ , TNF- α
130951MS	Q-Plex™ Mouse Cytokine-Screen IR (16-plex)	IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-10, IL-12p70, IL-17, MCP-1, IFN- γ , TNF- α , MIP-1 α , GM-CSF, RANTES

To place an order by phone/fax:

US Toll-free: 1 877-PBL-8881

International callers: +1 732-777-9123

OR

Fax a copy of your purchase order, including purchase order number or credit card information (MS/Visa only), contact name and telephone number to: +1 732-777-9141

To place an order by E-mail/Web:

Send an E-mail to sales@interferonsource.com with details of your purchase, your billing and shipping addresses (including Attn. to:), and your purchase order number or credit card information (MC/Visa only). You will receive an e-mail confirming your order. Alternatively, order via our convenient web form at www.interferonsource.com.

PBL InterferonSource, a division of Pestka Biomedical Laboratories, Inc. based in Piscataway, NJ, is the world's largest manufacturer and distributor of interferons and related products to the life sciences research market. Founded in 1990 by Dr. Sidney Pestka, PBL continues to deliver high quality products and timely services to provide researchers first-choice resources and solutions. Q-Plex and Quansys trademark under Quansys Biosciences, Utah, USA

