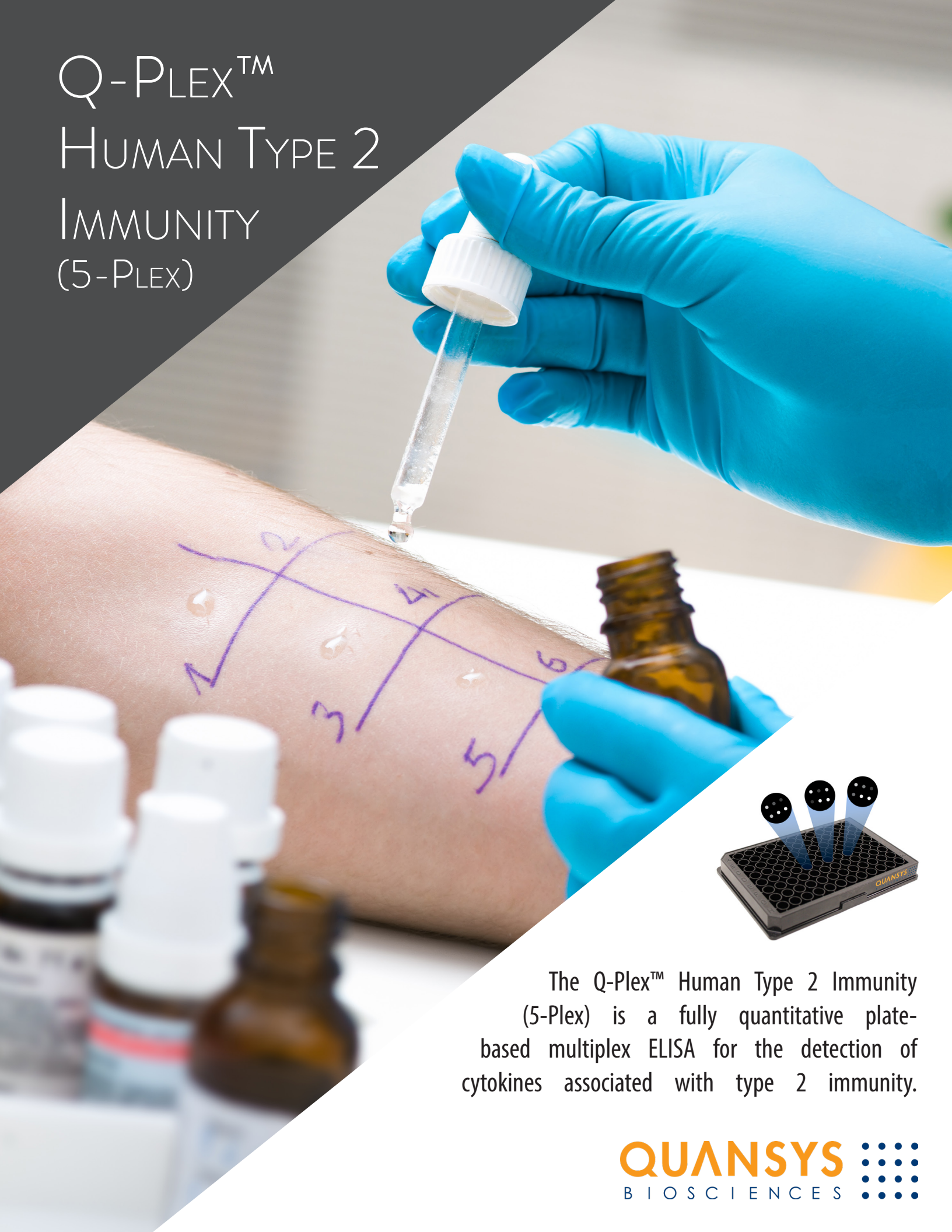
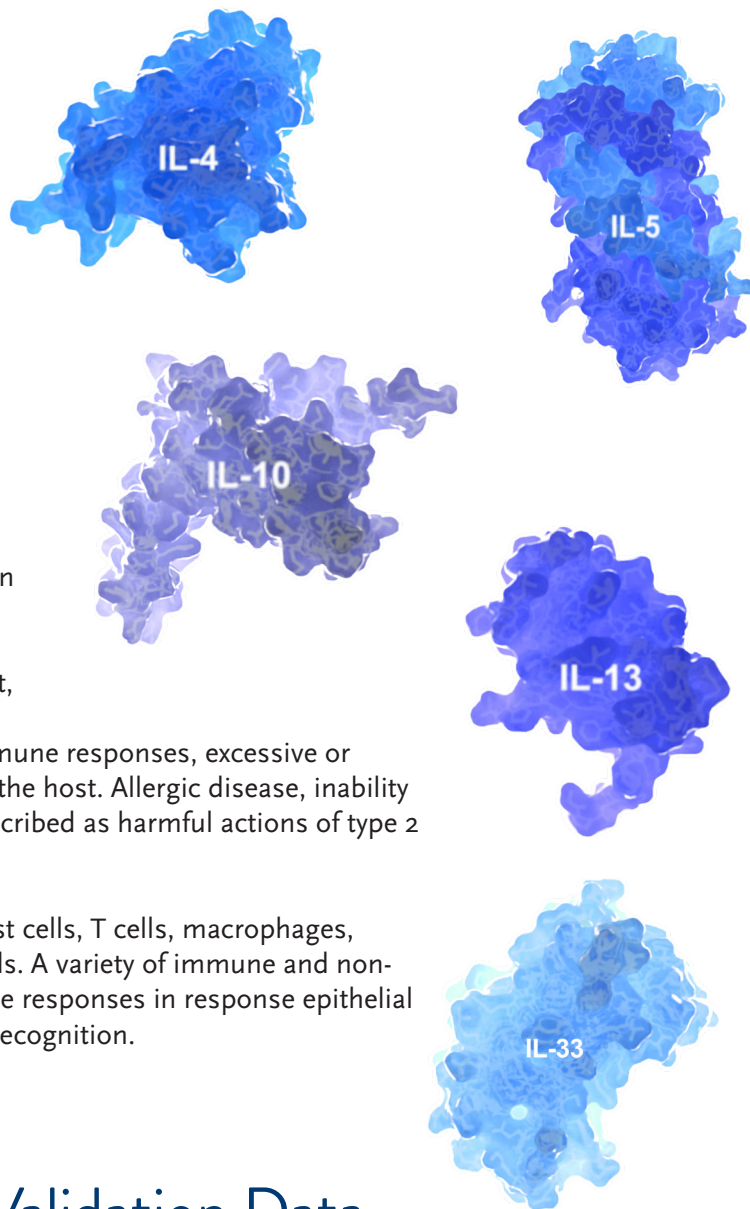


Q-PLEX™ HUMAN TYPE 2 IMMUNITY (5-PLEX)



The Q-Plex™ Human Type 2 Immunity (5-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with type 2 immunity.

Species Specificity	Human
Assay Type	Sandwich
Sample Type	Human Serum, EDTA/ Heparin Plasma
Assay Length	2.25 hours
Volume Required Per Well	Min. 25 µL
Detection Mention	Chemiluminescent
Multiplex Format	96-well solid plate
Catalog #	114233HU



Background

The Q-Plex™ Human Type 2 Immunity (5-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with type 2 immunity.

Functions of type 2 immunity include protecting the host, maintaining metabolic homeostasis, suppressing type 1, barrier defense, and tissue regeneration. Like other immune responses, excessive or chronic activation of type 2 immunity can be harmful to the host. Allergic disease, inability to defend against pathogens, and fibrosis have been described as harmful actions of type 2 immunity.

Producers of type 2 cytokines include macrophages, mast cells, T cells, macrophages, monocytes, basophils, eosinophils, and natural killer cells. A variety of immune and non-immune cells have been noted to produce type 2 cytokine responses in response epithelial damage, release of alarmins, and antigen presentation/recognition.

Example Performance and Validation Data

Analyte	Assay Type	Calibrator Range	Upper Limit of Quantification (ULOQ)	Lower Limit of Quantification (LLOQ)	Limit of Detection	Precision* (Inter-assay)	Precision* (Intra-assay)	Average* Linearity
IL-4	Sandwich	1,676 - 2.30 (pg/mL)	1,173 (pg/mL)	3.0 (pg/mL)	1.79 (pg/mL)	6%	4%	94%
IL-5	Sandwich	3,000 - 4.12 (pg/mL)	2,100 (pg/mL)	5.3 (pg/mL)	2.90 (pg/mL)	10%	5%	95%
IL-10	Sandwich	8,039 - 11.03 (pg/mL)	5,627 (pg/mL)	14.3 (pg/mL)	10.99 (pg/mL)	7%	4%	103%
IL-13	Sandwich	3,000 - 4.12 (pg/mL)	2,100 (pg/mL)	5.3 (pg/mL)	4.12 (pg/mL)	6%	5%	98%
IL-33	Sandwich	350 - 0.48 (pg/mL)	245 (pg/mL)	0.62 (pg/mL)	0.24 (pg/mL)	15%	6%	92%



Quansys Biosciences
365 North 600 West
Logan, UT 84321
www.quansysbio.com

Fax: 435-750-6869
Phone: 435-752-0531
Toll Free: 888-QUANSYS (782-6797)



Quansys is an ISO 9001:2015 and ISO 13485:2016 registered company and complies with GMP. Products are designed, developed, and manufactured according to the procedures outlined in our Quality Management System.

©2022 Quansys Biosciences Inc. All rights reserved. Q-View and Q-Plex are Trademarks of Quansys Biosciences Inc. ISO is a trademark of the International Standards Organization. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.