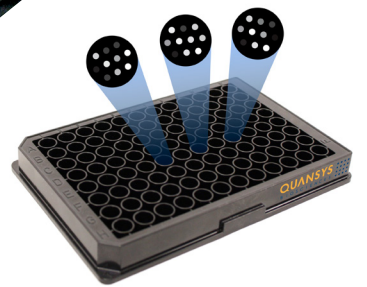


Q-PLEX™ HUMAN T CELL CYTOKINE HS (9-PLEX)



The Q-Plex™ Human T Cell Cytokine HS (9-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with T cell-mediated immunity.

Background

The Q-Plex™ Human T Cell Cytokine HS (9-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with T cell-mediated immunity.

T cells are one of the major lymphocyte arms of the adaptive immune system. Different subsets of T cells exist including Th1, Th2, Th17, and Tregs. Each subset of T cell is characterized by signature cytokine profiles as well as transcription factors and are specialized for protecting the host against different threats. T cell biology is studied in a variety of conditions including cancer, autoimmunity, allergy, and infectious diseases.

Species Specificity	Human
Assay Type	Sandwich
Sample Type	Human Serum, EDTA/ Heparin Plasma
Assay Length	4.5 hours
Volume Required Per Well	Min. 25 µL
Detection Mention	Chemiluminescent
Multiplex Format	96-well solid plate
Catalog #	111933HU

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
IFNγ	Sandwich	99 - 0.14 (pg/mL)	69.3 (pg/mL)	0.18 (pg/mL)	0.10 (pg/mL)	10%	3%	99%
IL-2	Sandwich	64 - 0.09 (pg/mL)	44.8 (pg/mL)	0.11 (pg/mL)	0.09 (pg/mL)	8%	6%	101%
IL-4	Sandwich	21 - 0.03 (pg/mL)	14.7 (pg/mL)	0.04 (pg/mL)	0.02 (pg/mL)	11%	4%	97%
IL-5	Sandwich	333 - 0.46 (pg/mL)	233 (pg/mL)	0.59 (pg/mL)	0.46 (pg/mL)	10%	7%	102%
IL-10	Sandwich	2,680 - 3.68 (pg/mL)	4.78 (pg/mL)	2.39 (pg/mL)	2.98 (pg/mL)	11%	3%	99%
IL-13	Sandwich	333 - 0.46 (pg/mL)	233 (pg/mL)	0.59 (pg/mL)	0.23 (pg/mL)	11%	5%	105%
IL-17a	Sandwich	5,589 - 7.67 (pg/mL)	3,912 (pg/mL)	9.97 (pg/mL)	6.21 (pg/mL)	9%	5%	104%
IL-21	Sandwich	250 - 0.34 (pg/mL)	175 (pg/mL)	1 (pg/mL)	0.17 (pg/mL)	12%	9%	95%
TNFα	Sandwich	661 - 0.91 (pg/mL)	462 (pg/mL)	1.18 (pg/mL)	0.46 (pg/mL)	12%	6%	100%



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.