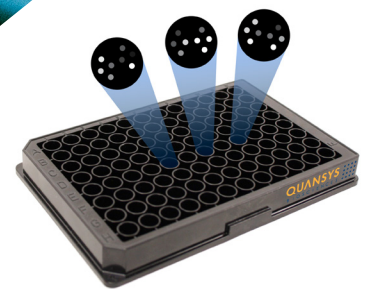


Q-PLEX™ HUMAN INNATE IMMUNITY (7-PLEX)



The Q-Plex™ Human Innate Immunity (7-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with innate immunity.

QUANSYS 
BIOSCIENCES

Background

The Q-Plex™ Human Innate Immunity (7-Plex) is a fully quantitative plate-based multiplex ELISA for the detection of cytokines associated with innate immunity.

The innate immune system includes anatomical barriers, chemical mediators such as the complement system, as well as immune cells capable of producing inflammatory cytokines. Innate immune cells are crucial to the host's defense against pathogens and for initiating immune responses.

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 Method	Chemiluminescent
Multiplex Format	96-well solid plate
Catalog #	111233HU

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
GMCSF	Sandwich	6,000 - 8.2 (pg/mL)	4,200 (pg/mL)	10.7 (pg/mL)	4.12 (pg/mL)	11%	7.20%	95%
IFNα	Sandwich	400 - 0.5 (pg/mL)	280 (pg/mL)	0.72 (pg/mL)	0.27 (pg/mL)	9%	5%	100%
IFNβ	Sandwich	19,000 - 26.1 (pg/mL)	21,850 (pg/mL)	16,150 (pg/mL)	18.57 (pg/mL)	11.7%	7.1%	110%
IFNγ	Sandwich	8,000 - 10.97 (pg/mL)	5,600 (pg/mL)	14.3 (pg/mL)	4.95 (pg/mL)	8%	6%	97%
IL-1α	Sandwich	4,000 - 5.49 (pg/mL)	2,800 (pg/mL)	7.1 (pg/mL)	5.43 (pg/mL)	6%	3%	86%
IL-1β	Sandwich	12,226 - 16.77 (pg/mL)	8,558 (pg/mL)	21.8 (pg/mL)	12.40 (pg/mL)	8%	6%	84%
IL-18	Sandwich	215 - 0.29 (pg/mL)	151 (pg/mL)	0.38 (pg/mL)	0.15 (pg/mL)	13%	4%	102%



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.