



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080235

Expiration Date: 31 OCT 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1310	pg/mL
IL-1 β	577	pg/mL
IL-2	1530	pg/mL
IL-4	228	pg/mL
IL-6	753	pg/mL
IL-8	580	pg/mL
IL-10	370	pg/mL
IL-12p70	518	pg/mL
IL-13	447	pg/mL
TNF- α	386	pg/mL

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 NOV 2019