



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080217

Expiration Date: 30 JUN 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1 330	pg/mL
IL-1 β	578	pg/mL
IL-2	1 550	pg/mL
IL-4	285	pg/mL
IL-6	752	pg/mL
IL-8	560	pg/mL
IL-10	383	pg/mL
IL-12p70	504	pg/mL
IL-13	500	pg/mL
TNF- α	342	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		25 JUL 2019