



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080348

Expiration Date: 30 NOV 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1500	pg/mL
IL-1 β	644	pg/mL
IL-2	1530	pg/mL
IL-4	258	pg/mL
IL-6	720	pg/mL
IL-8	601	pg/mL
IL-10	384	pg/mL
IL-12p70	488	pg/mL
IL-13	600	pg/mL
TNF- α	362	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	Nithya Santhanam	Quality		12 Jan 2023