



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080345

Expiration Date: 31 OCT 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1470	pg/mL
IL-1 β	620	pg/mL
IL-2	1450	pg/mL
IL-4	245	pg/mL
IL-6	710	pg/mL
IL-8	633	pg/mL
IL-10	394	pg/mL
IL-12p70	473	pg/mL
IL-13	566	pg/mL
TNF- α	364	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		04 JAN 2023