



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080216

Expiration Date: 30 JUN 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1280	pg/mL
IL-1 β	571	pg/mL
IL-2	1500	pg/mL
IL-6	755	pg/mL
IL-8	565	pg/mL
IL-10	373	pg/mL
IL-12p70	485	pg/mL
IL-13	486	pg/mL
TNF- α	320	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		23 DEC 2020