



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080342
Expiration Date: 30 SEP 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1700	pg/mL
IL-1 β	698	pg/mL
IL-2	1620	pg/mL
IL-4	248	pg/mL
IL-6	760	pg/mL
IL-8	674	pg/mL
IL-10	407	pg/mL
IL-12p70	498	pg/mL
IL-13	501	pg/mL
TNF- α	364	pg/mL

Reconstitution Volume: 1000 μ L

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		18 NOV 2022