



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080355
Expiration Date: 31 JAN 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1455	pg/mL
IL-1 β	463	pg/mL
IL-2	1480	pg/mL
IL-4	190	pg/mL
IL-6	760	pg/mL
IL-8	436	pg/mL
IL-10	368	pg/mL
IL-12p70	456	pg/mL
IL-13	545	pg/mL
TNF- α	290	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	Karen Hamilla	Quality		27 MAR 2023