



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080401
Expiration Date: 30 JUN 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	917	pg/mL
IL-1 β	1680	pg/mL
IL-2	2500	pg/mL
IL-4	1590	pg/mL
IL-5	889	pg/mL
IL-6	4650	pg/mL
KC/GRO	2050	pg/mL
IL-10	2790	pg/mL
IL-12p70	30700	pg/mL
TNF- α	591	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		26 JUL 2024