



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend

**Catalog Number:** C0049-2

**Lot Number:** A0080262

**Expiration Date:** 30 JUN 2025

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1560	pg/mL
IL-1 $\beta$	583	pg/mL
IL-2	1250	pg/mL
IL-4	213	pg/mL
IL-6	663	pg/mL
IL-8	575	pg/mL
IL-10	346	pg/mL
IL-12p70	447	pg/mL
IL-13	466	pg/mL
TNF- $\alpha$	338	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Carissa B. Hubbard	Quality		10 JUL 2020