



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Calibrator Blend

Catalog Number: C0084-2

Lot Number: A0080295

Expiration Date: 31 MAY 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A/F	6500	pg/mL
IL-17B	1130	pg/mL
IL-17C	2645	pg/mL
IL-17D	8645	pg/mL
IL-1RA	1095	pg/mL
IL-3	3260	pg/mL
IL-9	1575	pg/mL
TSLP	502	pg/mL

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 FEB 2022