



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Calibrator Blend

Catalog Number: C0084-2

Lot Number: A0080193

Expiration Date: 29 FEB 2024

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay							
	hu IL-17A/F	hu IL-17B	hu IL-17C	hu IL-17D	hu IL-1RA	hu IL-3	hu IL-9	hu TSLP
Reconstitution Calibrator (pg/mL)	5900	1210	2790	8780	905	2590	1620	494

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		06 MAR 2019