



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080177

Expiration Date: 31 DEC 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay								
	rt IFN- γ	rt IL-1 β	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- α
Reconstituted Calibrator (pg/mL)	5790	13300	1100	9230	13200	2770	27700	1610	1200

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	<i>Dawn Stollar</i>	08 JAN 2019