Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number:C0049-2Lot Number:A0080411Expiration Date:31 OCT 2029

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure	
IFN-γ	1410	pg/mL	
IL-1β	577	pg/mL	
IL-2	1410	pg/mL	
IL-4	220	pg/mL	
IL-6	751	pg/mL	
IL-8	603	pg/mL	
IL-10	385	pg/mL	
IL-12p70	507	pg/mL	
IL-13	534	pg/mL	
TNF-α	369	pg/mL	

Reconstitution Volume: 1000 µL

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	Laur Hamille	18 NOV 2024

COA-21479-A QA-FM-158-E