

EPX650-10065-901 ProcartaPlex Human Immune Monitoring Panel, 65-plex (As determined for):

(Assay Range determined for Standard Mix M Lot#144127000)

Analyte (Bead region)	Assay Range (pg/mL)
APRIL (88)	82.37–337400
BAFF (86)	8.35–34200
BLC (29)	12.23–50100
CD30 (84)	10.35–42400
CD40L (74)	4.76–19500
ENA-78 (82)	4.61–18900
Eotaxin (33)	1.65–6750
Eotaxin-2 (30)	4.39–18000
Eotaxin-3 (49)	2.23–9150
FGF-2 (75)	5.1–20900
Fractalkine (59)	2.47–10100
G-CSF (42)	9.86–40400
GM-CSF (44)	15.97–65400
Gro alpha (61)	3.1–12700
HGF (46)	3.44–14100
IFN-alpha (48)	4.69–19200
IFN-gamma (43)	13.13–53800
IL-1 alpha (62)	1.61–6600
IL-1 beta (18)	3.74–15300
IL-10 (28)	2.21–9050
IL-12p70 (34)	7.4–30300
IL-13 (35)	3.61–14800
IL-15 (65)	3.34–13700

Analyte (Bead region)	Assay Range (pg/mL)
IL-16 (70)	17.26–70700
IL-17A (36)	24.34–99700
IL-18 (66)	9.72–39800
IL-2 (19)	21.97–90000
IL-20 (81)	9.06–37100
IL-21 (72)	9.08–37200
IL-22 (76)	17.82–73000
IL-23 (63)	14.79–60600
IL-27 (14)	17.36–71100
IL-2R (9)	84.03–344200
IL-3 (73)	22.97–94100
IL-31 (37)	19.14–78400
IL-4 (20)	19.82–81200
IL-5 (21)	6.74–27600
IL-6 (25)	7.98–32700
IL-7 (26)	0.45–1850
IL-8 (27)	2.78–11400
IL-9 (52)	9.89–40500
IP-10 (22)	3.08–12600
I-TAC (57)	11.06–45300
LIF (15)	3.83–15700
MCP-1 (51)	3.83–15700
MCP-2 (8)	0.82–3350

Analyte (Bead region)	Assay Range (pg/mL)
MCP-3 (68)	4.83–19800
M-CSF (67)	65.45–268100
MDC (87)	13.43–55000
MIF (53)	0.78–3200
MIG (69)	8.64–35400
MIP-1 alpha (12)	3.05–12500
MIP-1 beta (47)	4.57–18700
MIP-3 alpha (56)	21.04–86200
MMP-1 (64)	5.13–21000
NGF beta (55)	5–20500
SCF (39)	3.39–13900
SDF-1 alpha (13)	56.86–232900
TNF alpha (45)	5.52–22600
TNF beta (54)	6.42–26300
TNF-R2 (85)	2.91–11900
TRAIL (58)	11.01–45100
TSLP (80)	4.71–19300
TWEAK (97)	709.72–2907000
VEGF-A (78)	6.84–28000