

Examples of Protein and Peptide Extracellular Mediators

ALCAM	Eot	GRO-g	IL-10	IL-21	MCP-4	b-NGF	TARC
ANG	Eot2	HB-EGF	IL-10 Rb	IL-3	M-CSF	NT-3	TGF-a
AR	Eot3	HCC-1	IL-11	IL-4	M-CSF R	NT-4	TGF-b1
BDNF	Fas	HCC-4	IL-12 p70	IL-5	MDC	OSM	TGF-b3
BLC	Fas Lig	HGF	IL-12 p40	IL-5 sRa	MIF	PARC	TIMP-1
VE-Cadherin	FGF-1	HVEM	IL-13	IL-6	MIG	PDGF Ra	TIMP-2
CCL28	FGF-2	I-309	IL-15	IL-6 sR	MIP-1a	PF4	TNF-a
CNTF	FGF-4	ICAM-1	IL-16	IL-7	MIP-1b	PIGF	TNF-b
CNTF Ra	FGF-6	ICAM-3	IL-17	IL-8	MIP-1d	Prolactin	TNF-RI
CT-1	FGF-7	IFN-a	IL-18	IL-9	MIP-3a	RANK	TNF-RII
CTACK	FGF-9	IFN-g	IL-1a	IP-10	MIP-3b	RANTES	TRAIL
CTLA-4	Flt-3 Lig	IFN-omega	IL-1b	I-TAC	MMP-1	SCF	TRAIL R1
CD27	Follistatin	IGF-I	IL-1ra	Leptin	MMP-10	SCF R	TRAIL R4
CD30	GCP-2	IGF-I R	IL-1 sRI	LIF	MMP-2	sCD23	uPAR
CD31	G-CSF	IGFBP-1	IL-1 sRII	sLIF R	MMP-7	SDF-1a	sVAP-1
CD40	GDNF	IGFBP-2	IL-2	Lymphotactin	MMP-9	SDF-1b	VEGF
DR6	GM-CSF	IGFBP-3	IL-2 sRa	MCP-1	MPIF-1	L-Selectin	VEGF R2
EGF	GRO-a	IGFBP-4	IL-2 sRb	MCP-2	MSP	sgp130	
ENA-78	GRO-b	IGF-II	IL-2 sRg	MCP-3	NAP-2	ST2	

