

# EXTRACELLULAR LIGANDS

www.signalchem.com

Cellular communication is influenced by both extra- and intra-cellular factors that can modulate or regulate specific signaling cascades in both positive and negative feedback mechanisms.

## CHEMOKINES

Product Name	Cat #
IP-10	I838-40N
SDF-1	S821-40N

## CYTOKINES

Product Name	Cat #
TNFB	T861-40N
TRAIL	T880-40N

## GROWTH FACTORS

Product Name	Cat #
BMP-2	B851-40N
BMP-4	B853-40N
IL-2	I862-40N

Product Name	Cat #
IL-8 (72AA)	I868-40N
Leptin	L833-40N
TGFB3	T833-40CN