





Human Receptor-Serine Kinases

STKR Family, Type1 Subfamily

ALK1/ACVRL1	
ALK2/ACVR1	
ALK4/ACVR1B	
ALK7/ACVR1C	
BMPR1A	
BMPR1B	
TGFβR1	

STKR Family, Type2 Subfamily

ACTR2/ACVR2A	
ACTR2B/ACVR2B	
BMPR2	
MISR2/AMHR2	
TGFβR2	