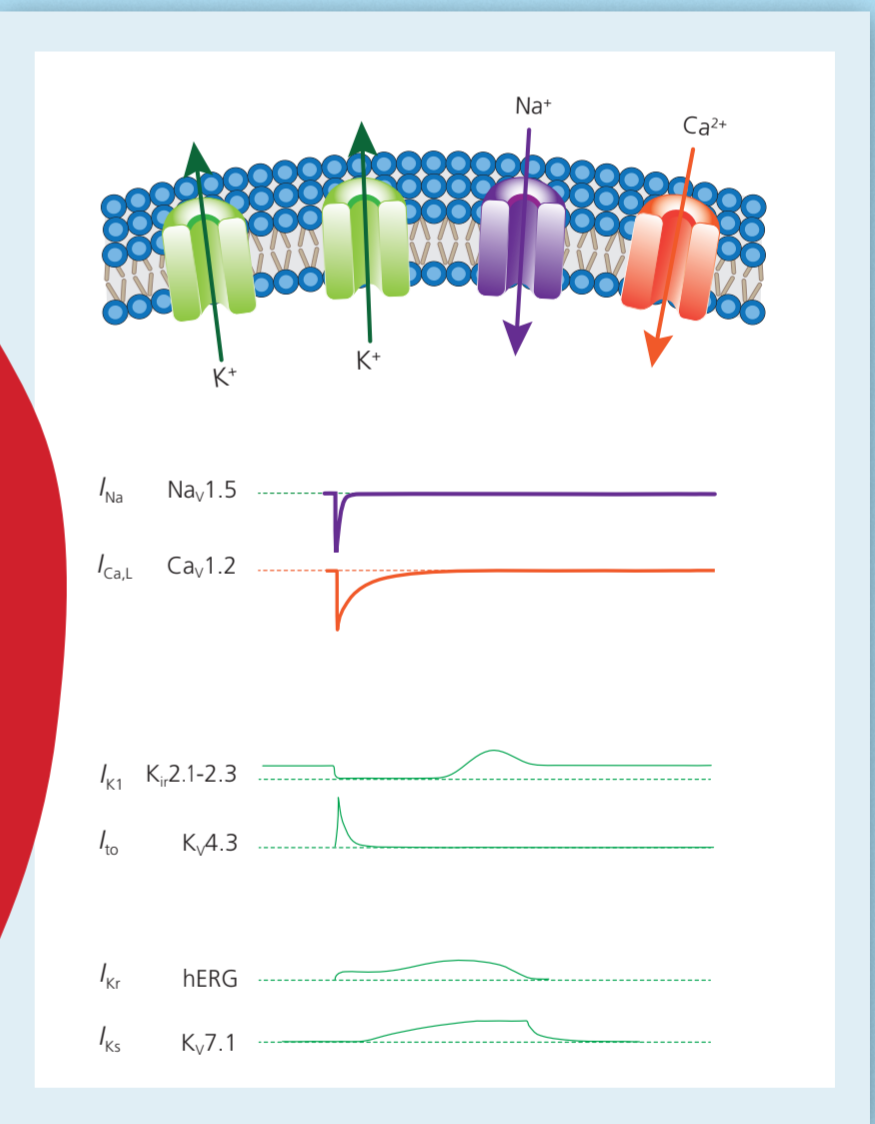
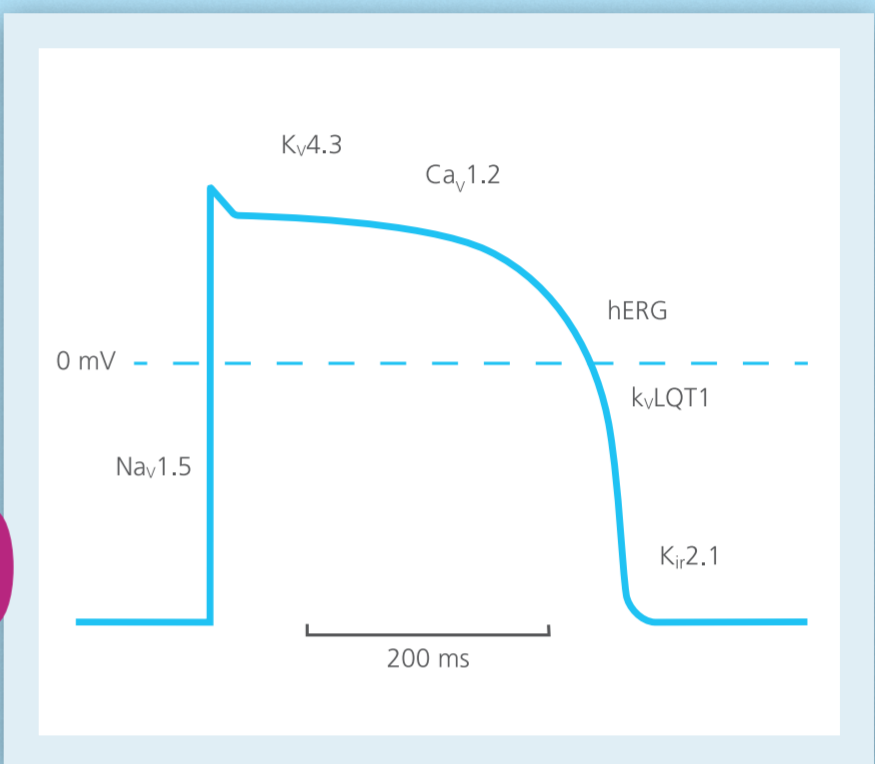
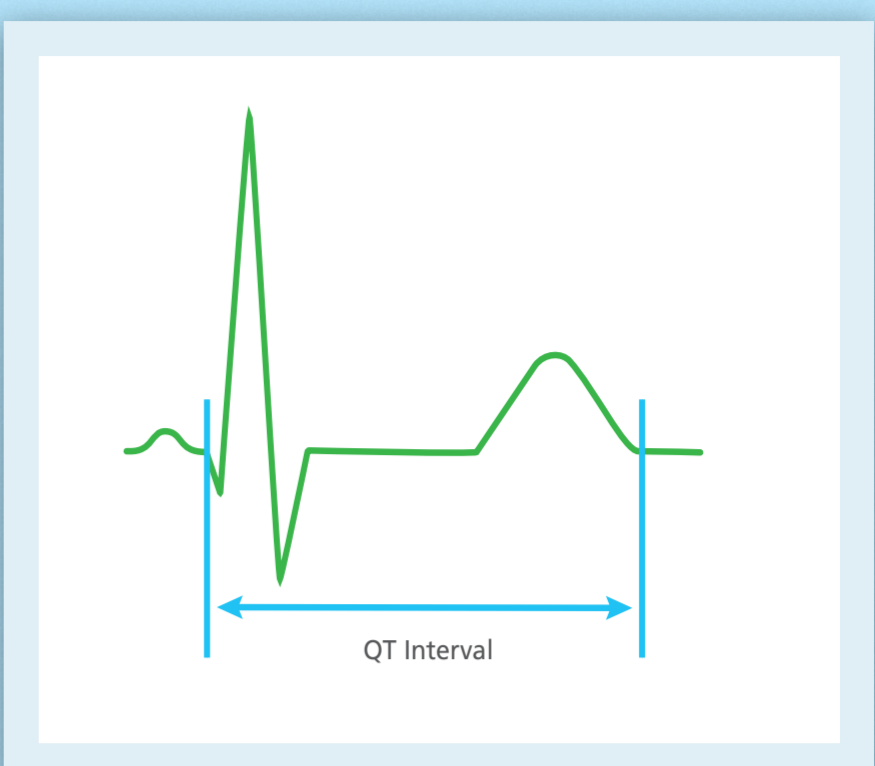
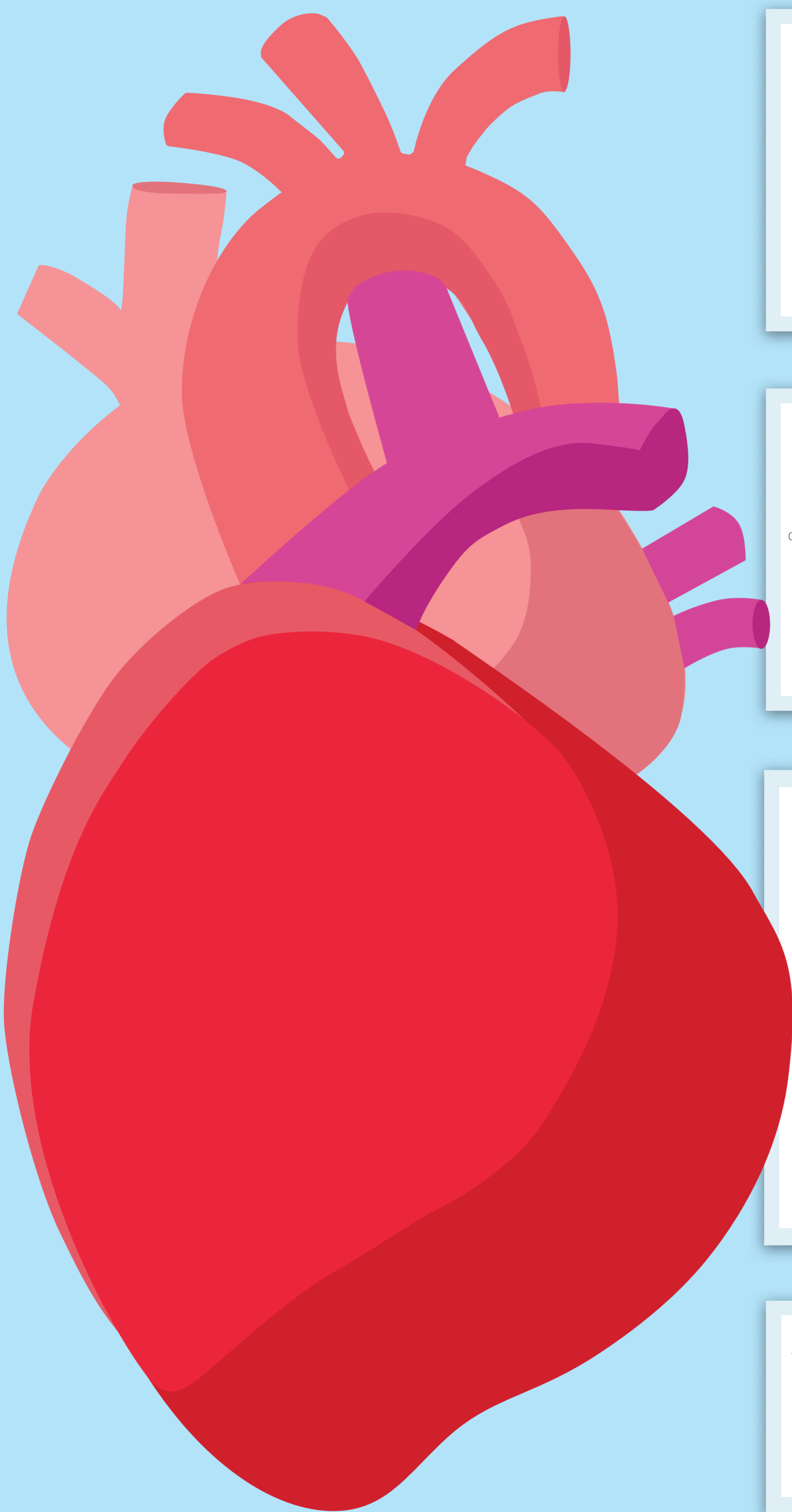


# CARDIAC ION CHANNELS



Ion Channel	Current	Gene
K <sub>v</sub> 11.1 (hERG)	$I_{Kr}$	KCNH2
Nav1.5	$I_{Na}$	SCN5A
Ca <sub>v</sub> 1.2	$I_{Ca,L}$	CACNA1C
K <sub>v</sub> 7.1 (KvLQT1)	$I_{Ks}$	KCNQ1
K <sub>v</sub> 4.3	$I_{to}$	KCND3
K <sub>i</sub> 2.1-2.3	$I_{K1}$	KCNJ2/12/4