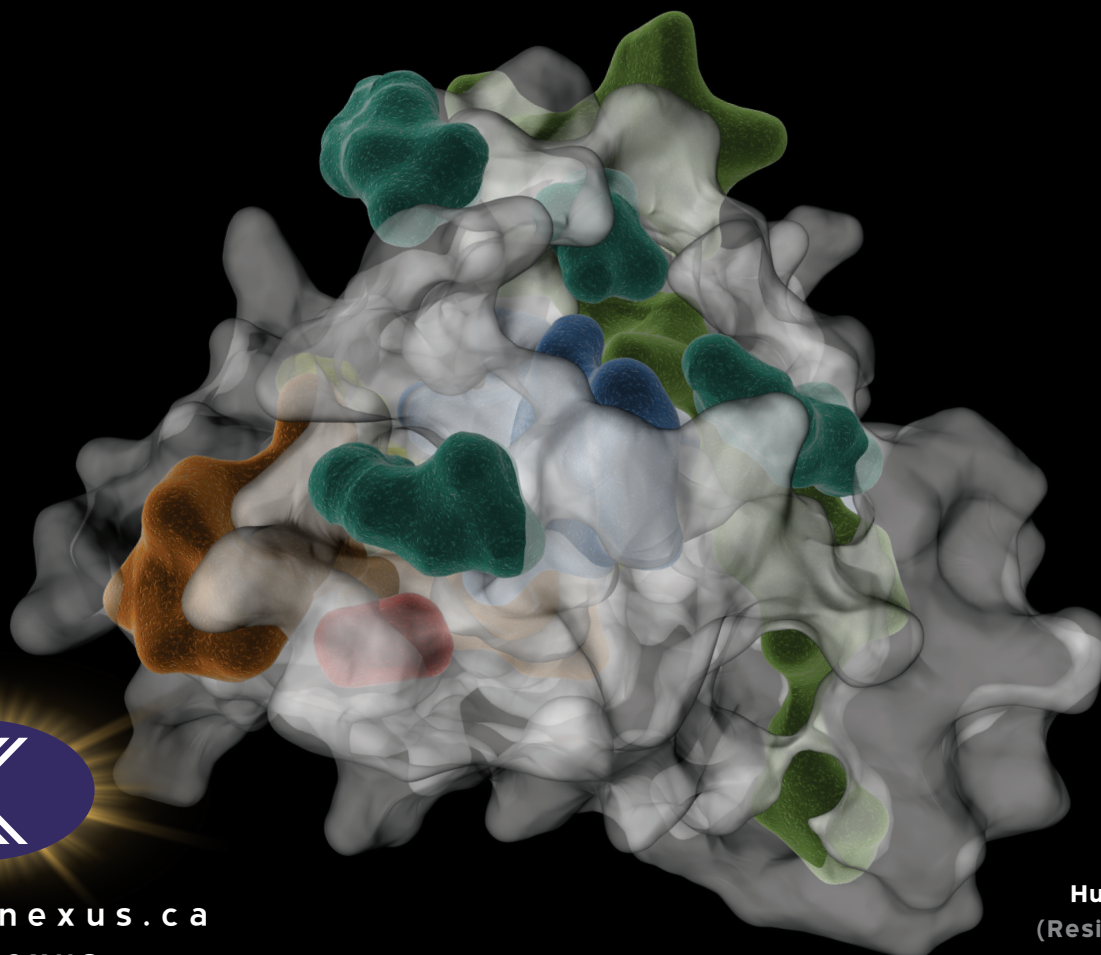


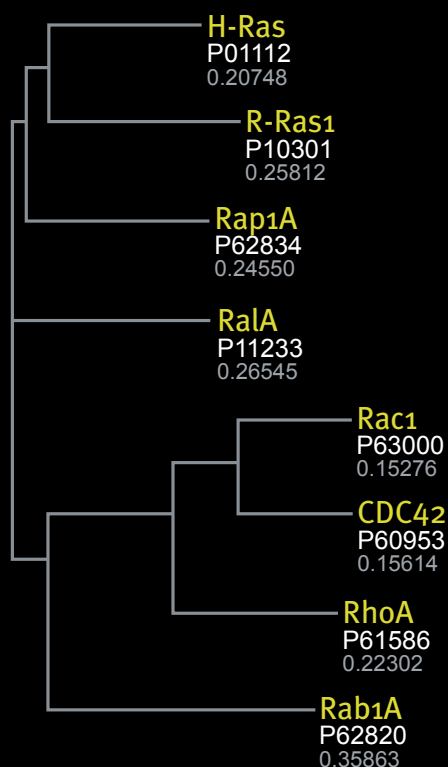


www.kinexus.ca
1-866-kinexus TOLL FREE PHONE



Human H-Ras
(Residues 1-166)
PDB: 121P

Ras-Related Small G Proteins



H-Ras	4	YKLVVVGAGGVGKSALT	TIQLIQNH	FVDEYDP	TIEDSYR	-KQVVIDGETCLLDILDTAGQEEYSAM
R-Ras1	30	HKLVVVGAGGVGKSALT	TIQFIQSY	FVSDYDP	TIEDSYT	-KICSDVGIPARLDILDTAGQEEFGAM
Rap1A	4	YKLVLGSGGVGKSALT	VQFVQGI	FVEKYDP	TIEDSYR	-KQVEVDCQCMLDILDTAGTEQFTAM
RalA	15	HKVIMVSGGVGKSALT	LQFMYDE	FVEDYEP	TKADSYR	-KKVVLGDGEEVQIDILDTAGQEDYAAI
Rab1A	12	FKLLLTIGDSGVGKSC	LLRFADDT	YTESYIS	TIGVDFK	IRTIELDGKTIKQLIWDTAGQERFRTI
Rac1	5	-KCVVVG	DGAVGKTCL	LISYTTNA	FPGEYIP	TVFDNYS-ANVMVDGKPVNLGLWDTAGQEDYDRL
CDC42	5	-KCVVVG	DGAVGKTCL	LISYTTNK	FPSEYVP	TVFDNYS-VTVMIGGEPYTLGLFDTAGQEDYDRL
RhoA	6	KKLVIV	GDGACGKTCL	LIVFSKDQ	FPEVYVP	TVFENYV-ADIEVDGKQVELALWDTAGQEDYDRL
H-Ras	68	RDQYMRTGEG	FLCVFA	INNTKSF	EDIHQYRE	QIKRVKSDDDVPMVLVGNKCDLAA-RTV
R-Ras1	94	REQYMRAGHG	FLLVFA	INDRQSF	NEVGKLF	TQILRVKDRDDFPVVLVGNKADLESQRQV
Rap1A	68	RDLYMKNQGQ	FALVYS	ITAQSTF	NDLQDLR	EQILRVKDTEDVPMILVGNKCDLEDERVV
RalA	79	RDNYFRSGEG	FLCVFS	ITEMESF	AATADFRE	QILRVKEDENVPFLLVGNKSDLEDKQV
Rab1A	77	TSSYYRGAHG	IIVVYD	VTDQESF	NNVQWLQ	EIDRYAS-ENVNKLVLVGNKCDLTTKKVV
Rac1	68	RPLSYPQTDV	FLICFSL	VSPA	SFENVR	AKWYPEVRHHC-PNTPILVGTKLDRDDKDT
CDC42	68	RPLSYPQTDV	FLVCFSL	VSPS	SFENVKEK	WVPEITHHC-PKTPFLLVGTQIDLRDDPST
RhoA	70	RPLSYPD	TDVILMCF	SIDSPD	SLENIPEK	WTPEVKHFC-PNVPIILVGNKCDLRNDEHT
H-Ras	126	ESRQAQDLARSYG	-IPYIET	SAKTRQ		150
R-Ras1	153	PRSEASAFGASHH	-VAYFEA	SAKLRL		177
Rap1A	127	GKEQGQNLAROWC	NAFL	ESAKSKI		152
RalA	138	SVEEAKNRAEQWN	-VNYVET	SAKTRA		162
Rab1A	87	DYTTAKEFADSLG	-IPFLET	SAKNAT		159
Rac1	126	IEKLKEKKLTPIT	YPQGLA	MAKEIGA		151
CDC42	126	IEKLAKNKQKPI	TETAEL	ARDLKA		151
RhoA	128	RRELAKMKQEPVK	PEEGRD	MANRIGA		153