

# Human Receptor Tyrosine Kinases - 1

## Alk Family

ALK

LTK



## Axl Family

TYRO3

AXL

MER/MERTK



## CCK4 Family

CCK4/PTK7



## DDR Family

DDR1

DDR2



## EGFR Family

EGFR

HER2/ErbB2

HER3/ErbB3

HER4/ErbB4



## Eph Family

EPHA1

EPHA2

EPHA3

EPHA4

EPHA5

EPHA6

EPHA7

EPHA8

EPHA10

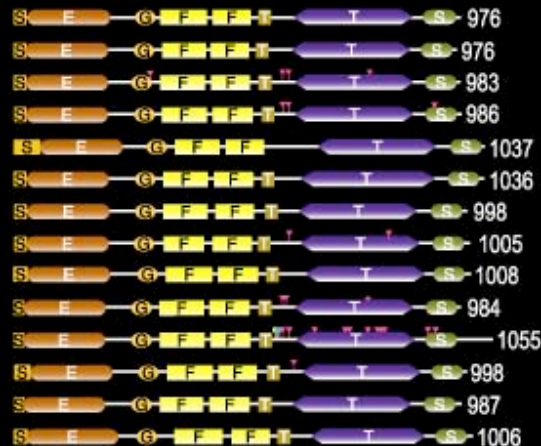
EPHB1

EPHB2

EPHB3

EPHB4

EPHB6



## FGFR Family

FGFR1

FGFR2

FGFR3

FGFR4



KINEXUS



www.kinexus.ca