

# Conservation of Human Phosphoproteome in Other Species

- The human and mouse phosphoproteomes are 46% different.
- Percent human phospho-sites conserved in mouse: 71.4
- Percent human phospho-sites conserved in frog: 14.4
- Percent human phospho-sites conserved in worm: 5.9
- Percent human phospho-sites conserved in yeast: 1.7