

Human Phosphoproteome Number of Phospho-Site Connections

- Likely number of human phosphosites: ~1,000,000
- Number of proteins: ~30,000
- Average number of sites per protein: ~33
- Average number of amino acids per protein: 500
- Frequency of phosphorylation sites: 1 every 15 aa
- Average number phospho-sites per kinase: ~2000
- Average number phospho-sites per phosphatase: ~7250
- Minimum number of connections in phosphorylation networks: ~2,000,000
- Likely number of connections in phosphorylation networks: ~10,000,000