

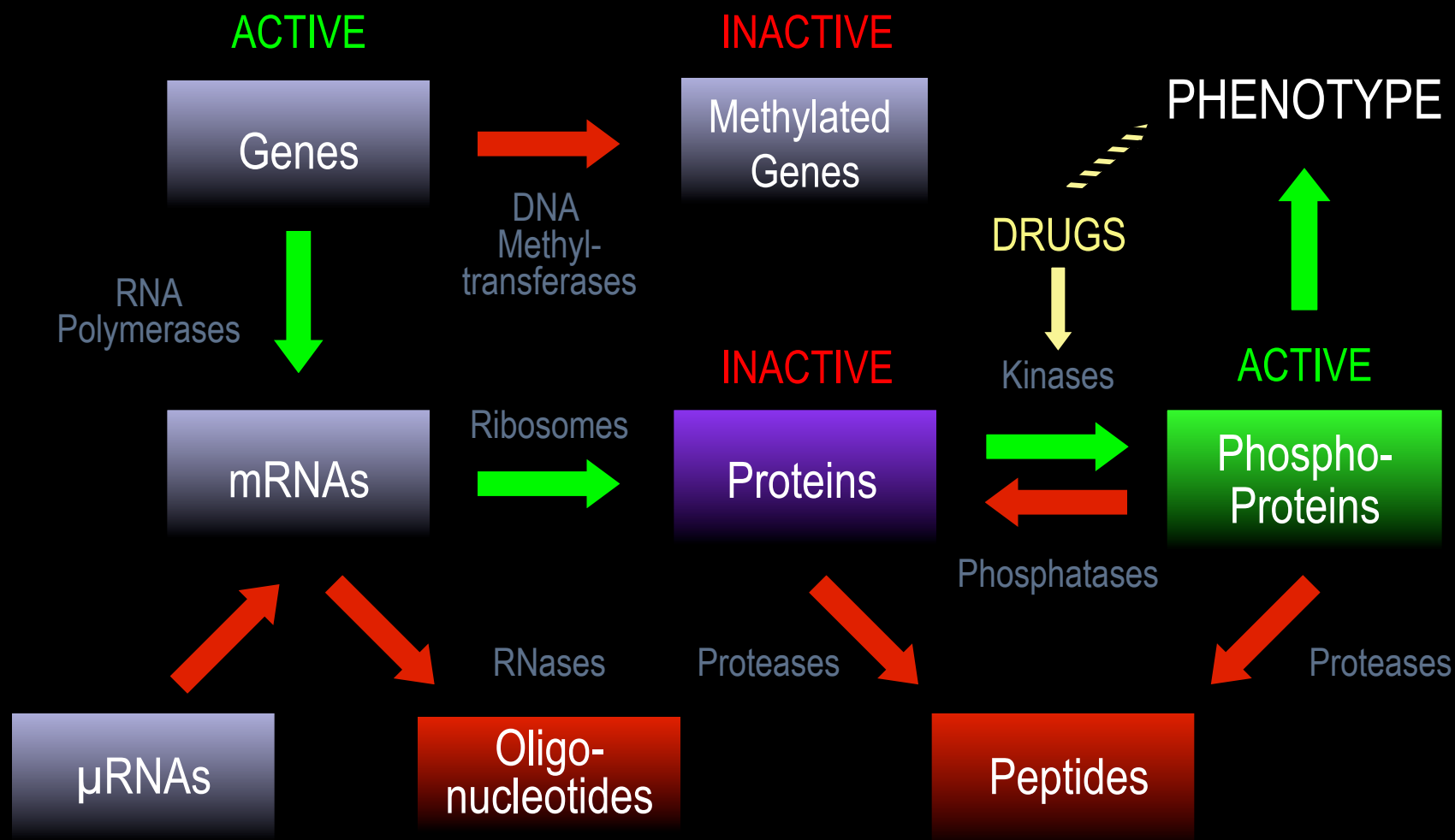
# Phosphoproteomics and the Origin and Operations of the Kineome

**Presented by Steven Pelech, Ph.D.**

Professor, Department of Medicine,  
University of British Columbia  
President & CSO, Kinexus Bioinformatics Corp.



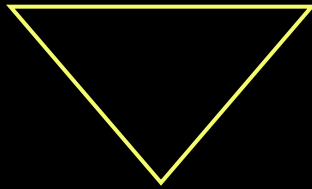
Presented at the OMICS Group 4<sup>th</sup> International Conference on  
Proteomics & Bioinformatics in Northbrook, IL, USA, 2014 Aug. 5



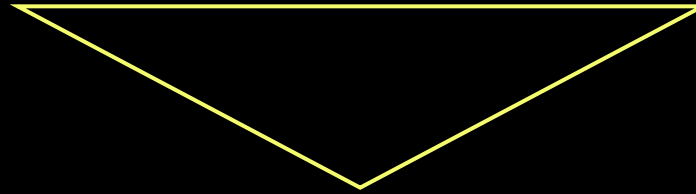
Common Name	Scientific Name	#Genes	#Protein Kinases
Baker's yeast	<i>Sacchar. cerevisiae</i>	6,300	122
Cat	<i>Felis silvestris</i>	20,285	
Chicken	<i>Gallus gallus</i>	23,000	546
Chimpanzee	<i>Pan troglodytes</i>	30,000	587
Cow	<i>Bos taurus</i>	22,000	
Dog	<i>Lupus familis</i>	40,000	656
Fruit fly	<i>Dros. melanogaster</i>	14,000	319 (232)
Honey bee	<i>Apis mellifera</i>	10,157	
Human	<i>Homo sapiens</i>	21,500	>536 (505)
Maize	<i>Zea mays</i>	32,000	
Mouse	<i>Mus musculus</i>	23,800	510 (527)
Nematode worm	<i>Caenorhab. elegans</i>	22,900	437 (434)
Poplar tree	<i>Populus trichocarpa</i>	45,555	
Rat	<i>Rattus norvegicus</i>	30,000	521
Red bread mold	<i>Neurospora crassa</i>	10,000	103
Rhesus Macaque	<i>Macaca mulatta</i>	18,296	821
Rice	<i>Oryza sativa</i>	56,000	1429
Sea urchin	<i>Strongylocentrotus purpuratus</i>	23,300	353
Thale Cress	<i>Arabidopsis thaliana</i>	27,000	1049
Tiger blowfish	<i>Takifugu rubripes</i>	31,000	519
Zebra Danio	<i>Brachydanio rerio</i>	24,200	



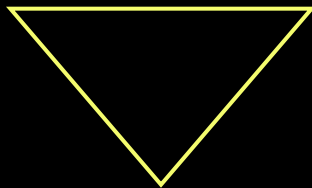
PROTEOME



~21,500 total protein genes

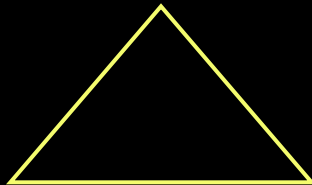


KINEOME



>536 protein kinase genes  
>1000 catalytically active protein  
kinase isoforms

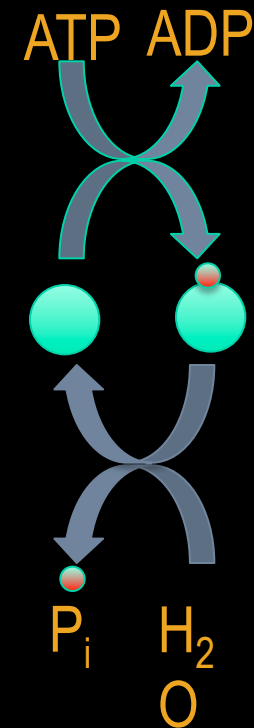
PHOSPHO-  
PROTEOME



~1,000,000 predicted  
phosphorylation sites (>180,000  
experimentally confirmed)

PHOSPHATOME

>156 protein phosphatase genes



## Statistics on the Human Phosphoproteome

- Number of human phosphosites known: >180,000
- Number of human phosphoproteins identified: >19,800
- Percent serine phosphorylation sites: 53.5%\*
- Percent threonine phosphorylation sites: 25.2%\*
- Percent tyrosine phosphorylation sites: 21.3%\*

\*Based on 159,540 human phosphosites identified by mass spectrometry



Based on 60,000 human phosphosites identified by mass spectrometry without specific enrichment of tyrosine phosphorylation

- Percent serine phosphorylation sites: 76%
- Percent threonine phosphorylation sites: 20%
- Percent tyrosine phosphorylation sites: 4%
- Number of tyrosine phosphosites known: >36,194

Minimum number of human phosphosites:  $(36,000 \times 25 =)$  900,000

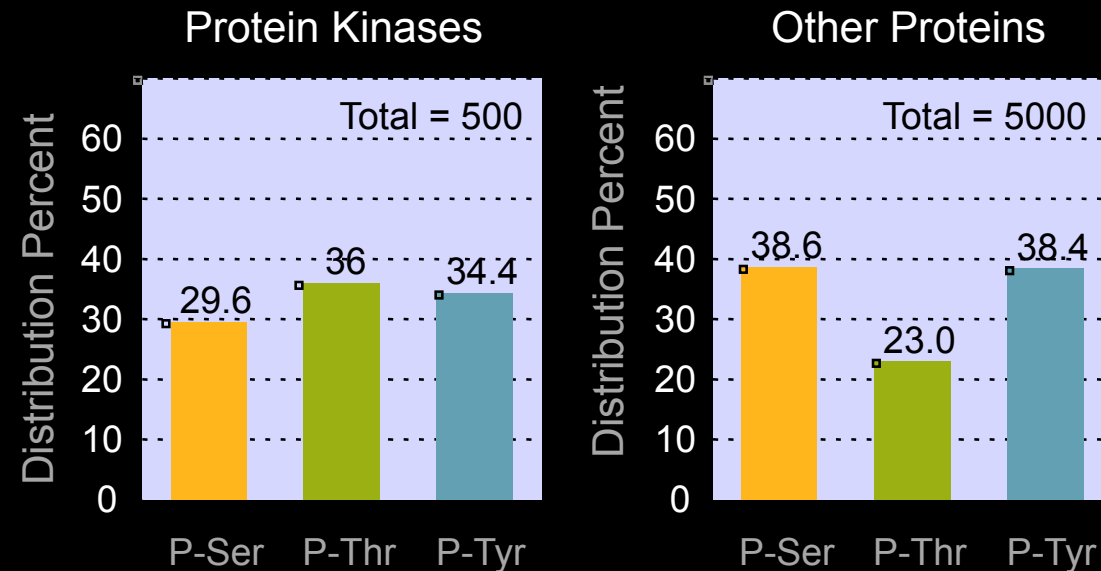
Probable number of human phosphosites: 1,000,000



	Selected Amino Acid		
	Serine	Threonine	Tyrosine
Percentage of total phosphosites	76%	20%	4%
Estimated number of total phosphosites (assume 1,000,000 total sites)	760,000	200,000	40,000
Estimated number of phosphosites per average protein (assume 21,500 proteins encoded by human genome)	35.3	9.30	1.86
Phosphosite frequency - Average length of peptide in amino acids with a phosphosite	19.9	75.6	378
Frequency of selected amino acid in general in human proteins	7.1%	6.0%	3.2%
Number of selected amino acids per average protein (average protein is 703)	49.91	42.18	22.50
Percent phosphosites per total available selected amino acids	70.7%	22%	8.3%
Ratio phosphosites per total available selected amino acids	2/3	1/5	1/12



### Most Evolutionary Conserved Human Phosphosites

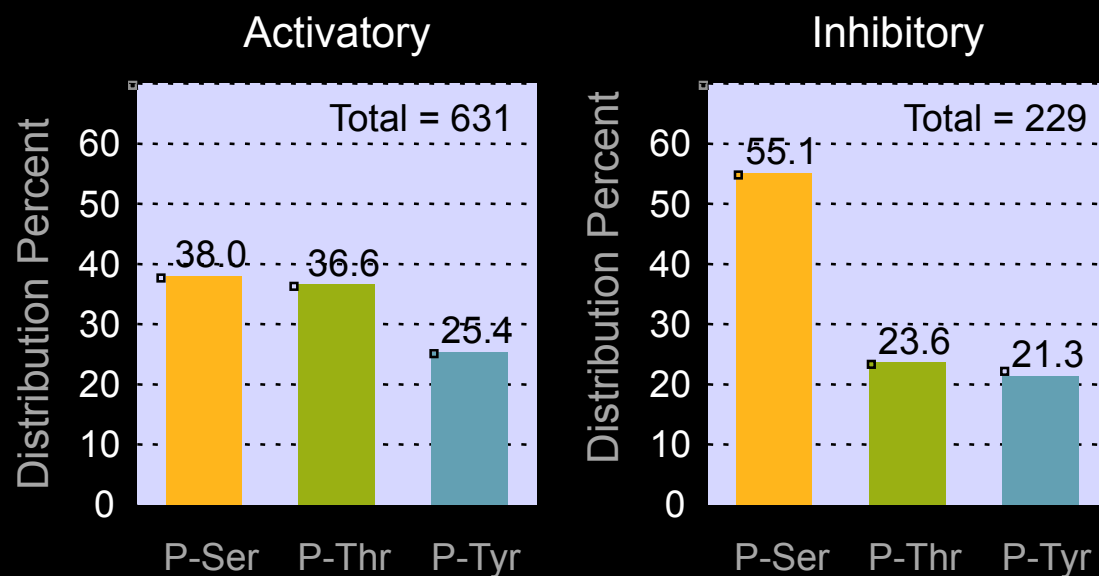


- About 9% of the most conserved phosphosites are found in protein kinases, which represent about 2.5% of all of the human protein genes
- While about 76% of phosphosites in general are P-Ser, in the most conserved phosphosites in protein kinases, P-Thr is the most prevalent





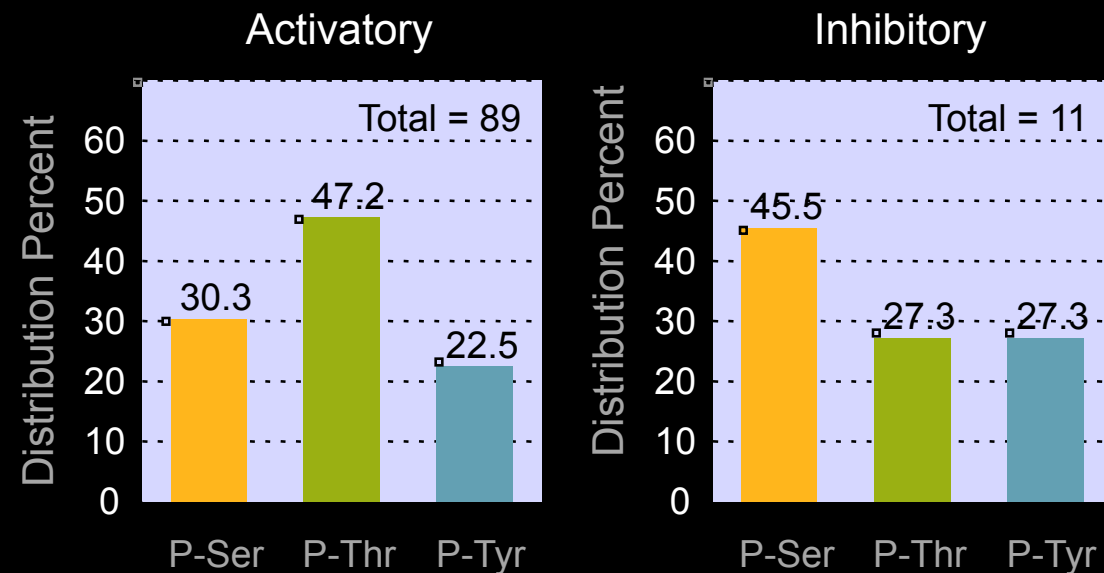
### Functional Human Kinase Phosphosites



- In general, functional phosphosites are more likely to be activatory than inhibitory – 2.8:1 ratio



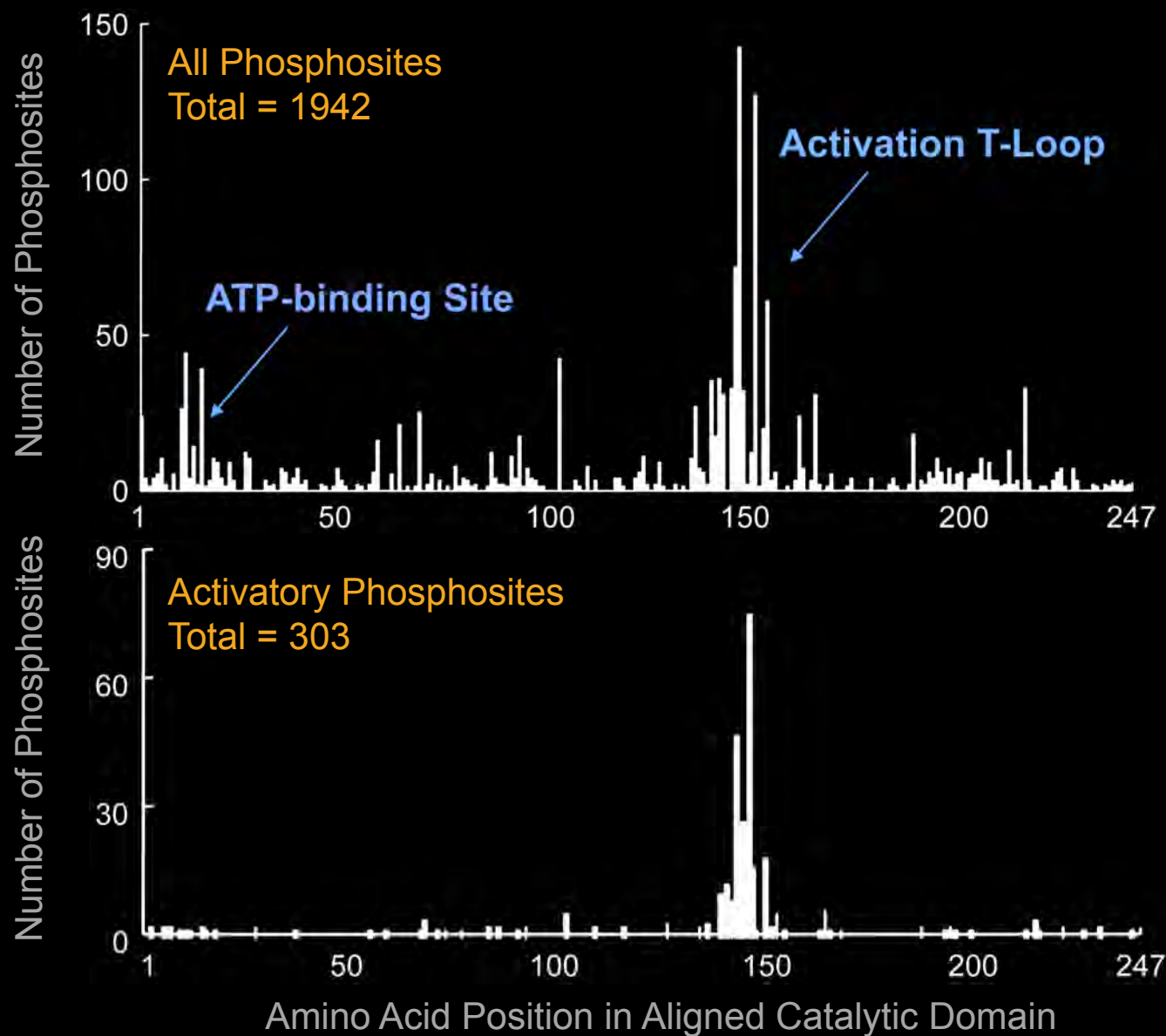
### Most Conserved Functional Human Kinase Phosphosites



- Functional phosphosites that are conserved are more likely to be activatory than inhibitory in kinases – 8:1 ratio



### Distribution of Phosphosites





## Phosphosites Between Kinase Subdomains VII and VIII

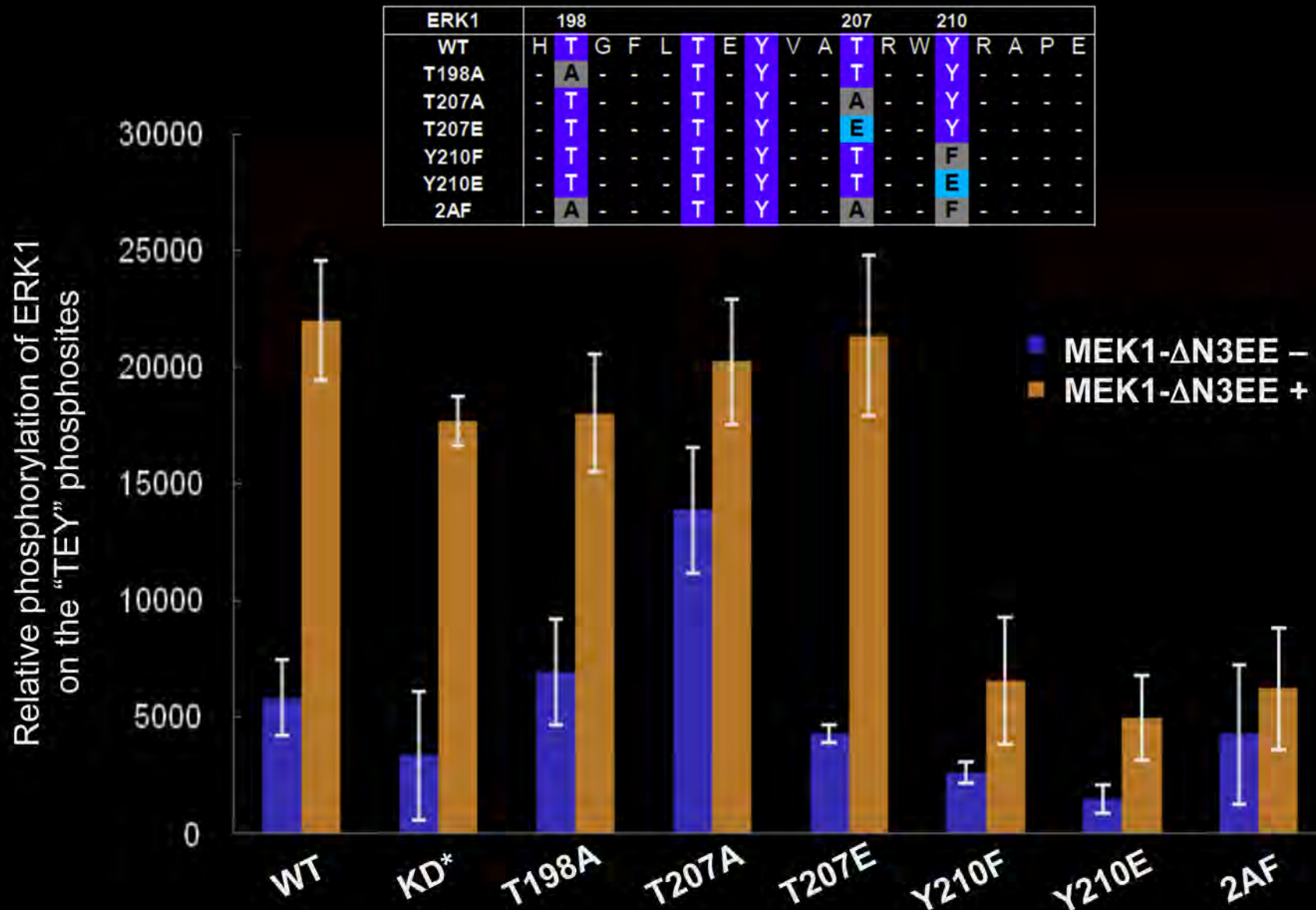
	144	145	146	147	148	149	150	151	152	153	154	155	156	157	158	159	160	161
CDK1	P	I	R	T	Y	T	H	E	V	V	T	L	W	Y	R	S	P	E
CDK2	P	V	R	T	Y	T	H	E	V	V	T	L	W	Y	R	A	P	E
GSK3	R	G	E	P	N	V	S	Y	I	C	S	R	Y	Y	R	A	P	E
ERK1	H	T	G	F	L	T	E	Y	V	A	T	R	W	Y	R	A	P	E
p38 alpha	T	D	D	E	M	T	G	Y	V	A	T	R	W	Y	R	A	P	E
JNK1	T	S	F	M	M	T	P	Y	V	V	T	R	Y	Y	R	A	P	E
PKC alpha	D	G	V	T	T	R	T	F	C	G	T	P	D	Y	I	A	P	E
PKC delta	G	E	S	R	A	S	T	F	C	G	T	P	D	Y	I	A	P	E
RSK1	H	E	K	K	A	Y	S	F	C	G	T	I	E	Y	M	A	P	E
MAPKAPK2	S	H	N	S	L	T	T	P	C	Y	T	P	Y	Y	V	A	P	E
CAMK1	P	G	S	V	L	S	T	A	C	G	T	P	G	Y	V	A	P	E
PIM1	K	D	T	V	Y	T	D	F	D	G	T	R	V	Y	S	P	P	E
AKT2	D	G	A	T	M	K	T	F	C	G	T	P	E	Y	L	A	P	E
PLK1	D	G	E	R	K	K	T	L	C	G	T	P	N	Y	I	A	P	E

Purple, confirmed activation sites;  
 Blue, confirmed phosphosites with unknown function;  
 Grey, potential phosphosites based on similarities

- Kinase Catalytic Domain residues 149-151 most commonly feature activatory phosphosites



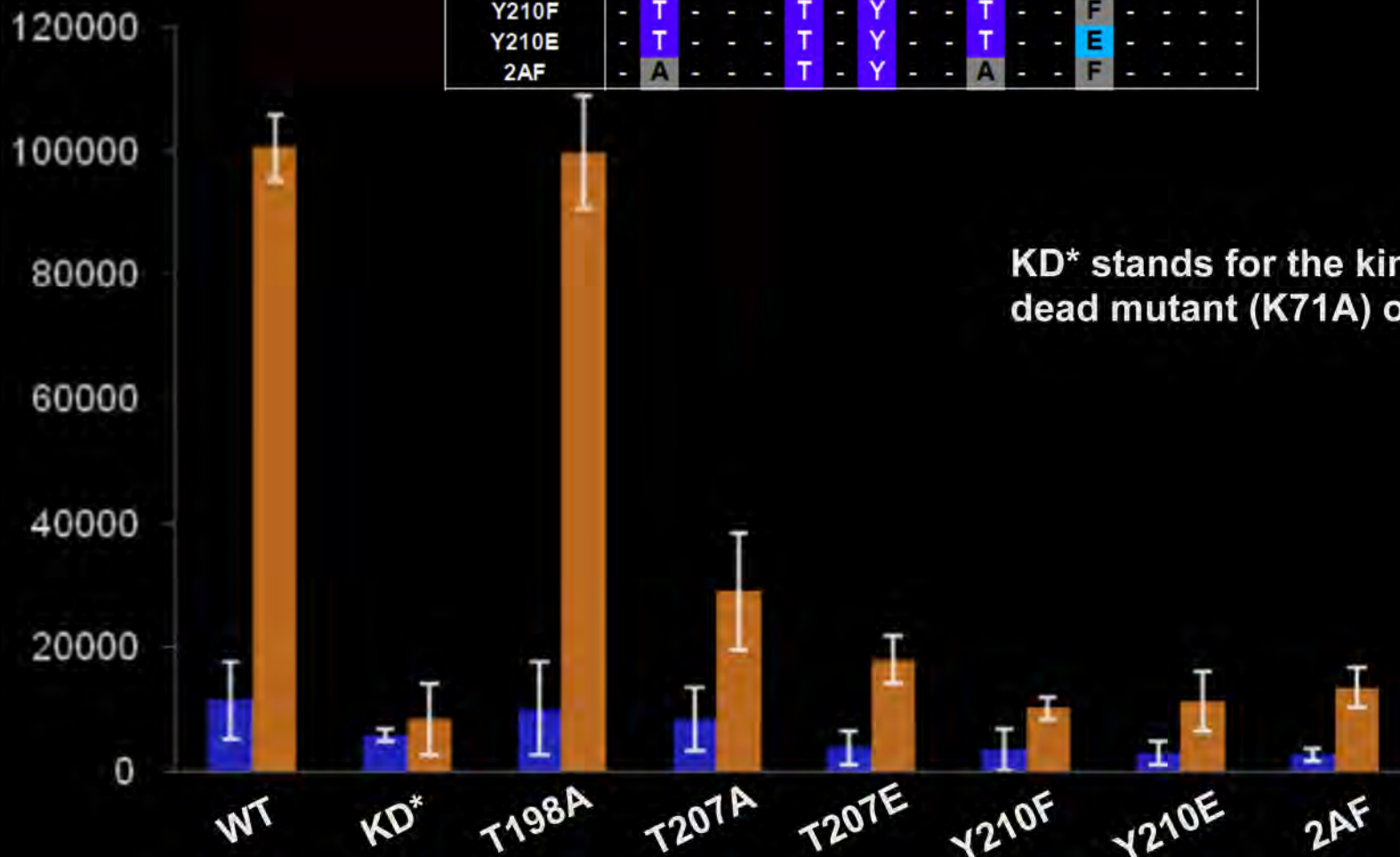
### Role Phosphosites Between Kinase Subdomains VII and VIII





### Phosphorylation of Myelin Basic Protein by ERK1

ERK1	198								207				210					
WT	H	T	G	F	L	T	E	Y	V	A	T	R	W	Y	R	A	P	E
T198A	-	A	-	-	-	T	-	Y	-	-	T	-	-	Y	-	-	-	-
T207A	-	T	-	-	-	T	-	Y	-	-	A	-	-	Y	-	-	-	-
T207E	-	T	-	-	-	T	-	Y	-	-	E	-	-	Y	-	-	-	-
Y210F	-	T	-	-	-	T	-	Y	-	-	T	-	-	F	-	-	-	-
Y210E	-	T	-	-	-	T	-	Y	-	-	T	-	-	E	-	-	-	-
2AF	-	A	-	-	-	T	-	Y	-	-	A	-	-	F	-	-	-	-



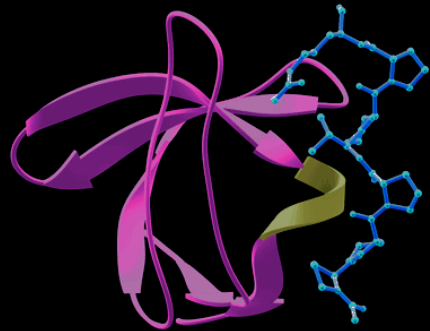


## Amino Acids Commonly Surrounding Serine Phosphosites

	-7	-6	-5	-4	-3	-2	-1	+1	+2	+3	+4	+5	+6	+7		
Ser	15	15	6	43	3	55	18	-44	40	20	36	17	20	15	S	Phospho-
Pro	6	-1	4	-6	-18	17	-1	326	8	4	-15	9	-8	6	P	serine
Glu	-4	-1	-13	5	-22	-20	-33	-20	40	68	20	20	11	16	E	at 0 position
Arg	14	23	50	20	179	65	-1	-57	-23	-3	-18	6	11	-5	R	
Ala	14	12	6	-8	-4	15	16	-42	1	8	-6	5	4	4	A	Based on
Glu	0	-5	-8	-5	-2	-22	27	-36	2	2	-7	-15	-3	-9	G	4035 human
Leu	-15	-6	26	-10	-10	-19	31	-18	-23	-16	4	-5	2	3	L	sites
Asp	-12	-21	-16	-8	-13	-15	26	32	25	18	5	16	6	2	D	
Lys	24	21	0	12	12	-24	-27	-64	-15	3	-4	-4	14	2	K	% Change
Thr	-11	-3	-21	-7	-21	6	-21	-55	-1	-27	5	-9	-6	-14	T	from
Val	-17	-12	-18	-21	-27	-32	-14	-23	13	-30	-9	-5	-10	-3	V	Expected
Gln	19	7	3	8	-4	-3	-22	-5	-7	-21	-5	-4	0	-4	Q	
Asn	-18	-11	-28	-16	-22	-18	8	-46	-29	-6	-14	-22	-16	-17	N	>100
Ile	-8	-14	7	-25	-32	-24	-7	-42	-30	-40	10	-18	-23	-5	I	50 to 100
Phe	-25	-13	-23	-4	-29	-43	1	12	-58	-33	-14	-14	-27	-23	F	25 to 49
Tyr	-24	-28	-27	-27	-38	-22	-27	-43	-34	-28	-33	-20	-14	-12	Y	0 to 24
His	-15	15	13	-24	-26	-40	-8	-46	-47	-27	-16	-2	-15	-4	H	-25 to -1
Met	-22	-12	15	-25	-45	-36	-3	-28	-37	-24	0	-15	-21	-7	M	-50 to -26
Cys	-4	-28	-32	-11	-28	5	-7	-63	-14	-25	-25	-34	-29	-21	C	-100 to -51
Trp	3	-52	-27	-42	-49	-49	-49	-42	-45	-37	-15	-44	-40	9	W	

Percentage Changes from Expected Random Frequency

- Proline appears at the +1 position in ~26% of P-Ser and P-Thr sites



**SH3 Domain**

e.g. **SEM5**

**Left-hand  
poly-Pro type II helix**

**P-X-X-P**

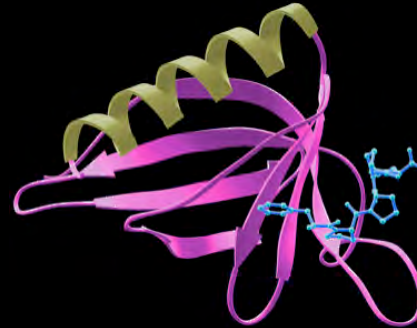


**WW Domain**

**Pin1**

**Pro-Rich**

**Pho-S/Pho-T  
P-P-X**



**EVH1 Domain**

**EVL**

**Pro-Rich**

**E/D-F-P-P-P-P-X-D/E**



**GYF Domain**

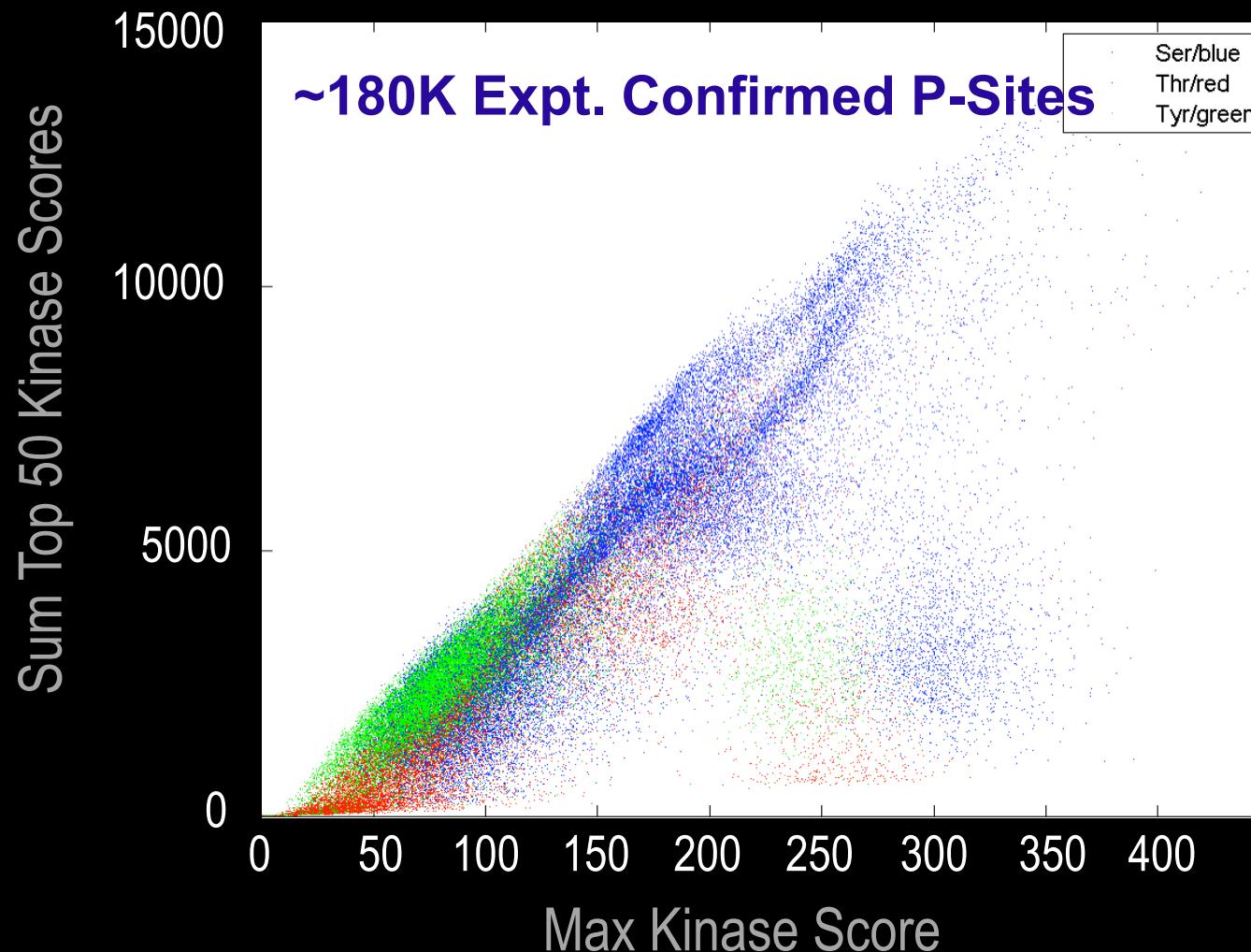
**CDC2BP2**

**Pro-Rich**

**P-P-P-P-G-H-R**

Poly-proline-rich binding domains mediate protein-protein interactions.  
Phosphorylation may regulate these interactions.





Very few phosphorylation sites appear to be targeted with high specificity by only a few kinases. Most phosphorylations are not functionally important.

- PEST sequences are rich in Pro, Glu, Ser and Thr
- PEST sequences are sites of known protein degradation  
e.g. cyclins C, D and E; PKC isoforms
- ~ 2,732 PEST sequences of 15 aa length with 3 or more prolines identified in human proteome (~18,619 sequences of 15 aa length with 2 or less prolines also identified - known as ST-rich sites)
- ~42% of PEST sequences appear to be phosphorylatable, compared to ~26% of ST-rich sites, despite more S and T in ST-rich sites
- Phosphorylation may serve to activate PEST sequences  
e.g. plant FRQ and CK1 phosphorylation

## Relationship between Protein Phosphorylation and Mass

Based on 80 phospho-sites using immunoblotting (data from [www.kinet.ca](http://www.kinet.ca))

### ■ Serine Phosphosites

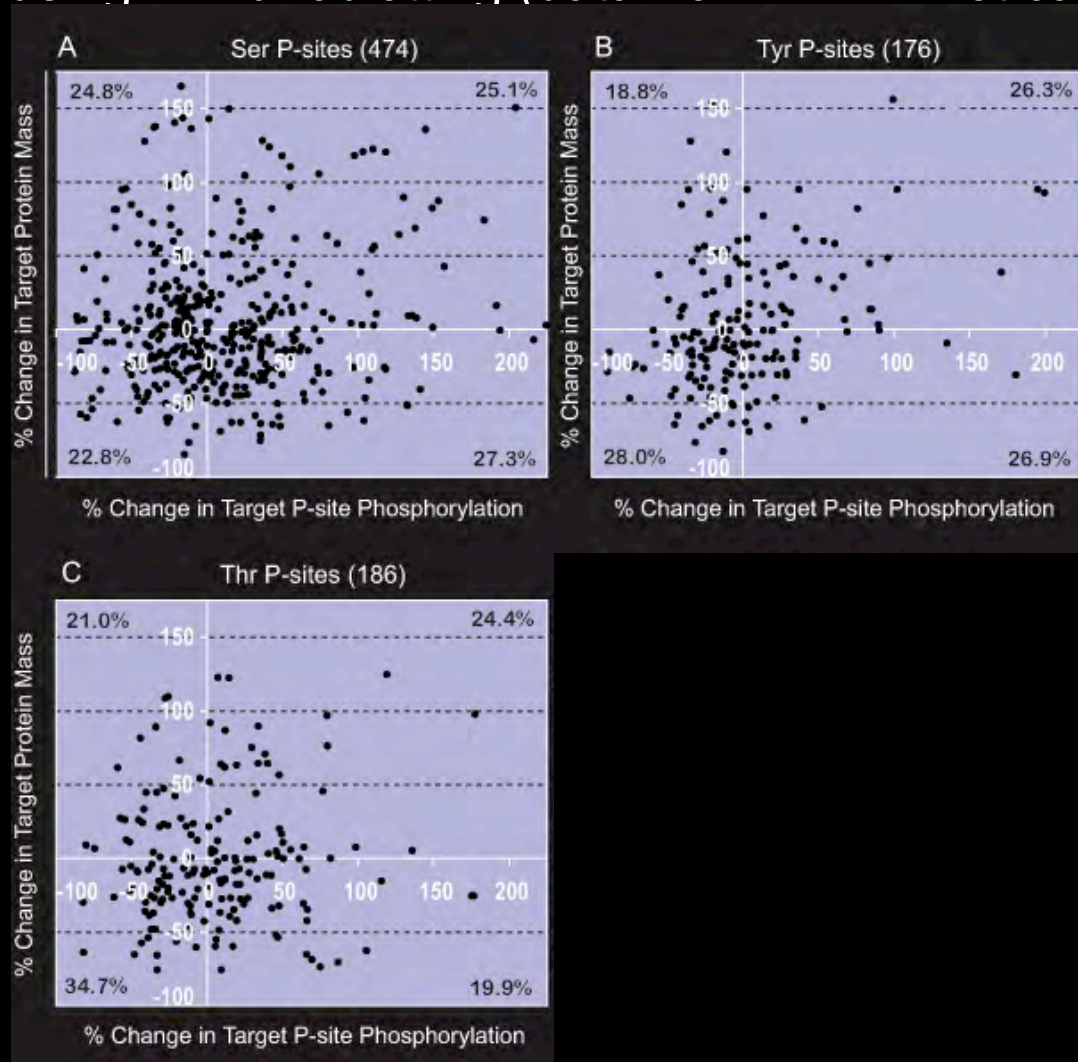
- Positive correlation = 47.9 %
- Negative correlation = 52.1 %

### ■ Threonine Phosphosites

- Positive correlation = 59.1 %
- Negative correlation = 40.9 %

### ■ Tyrosine Phosphosites

- Positive correlation = 45.7 %
- Negative correlation = 54.3 %

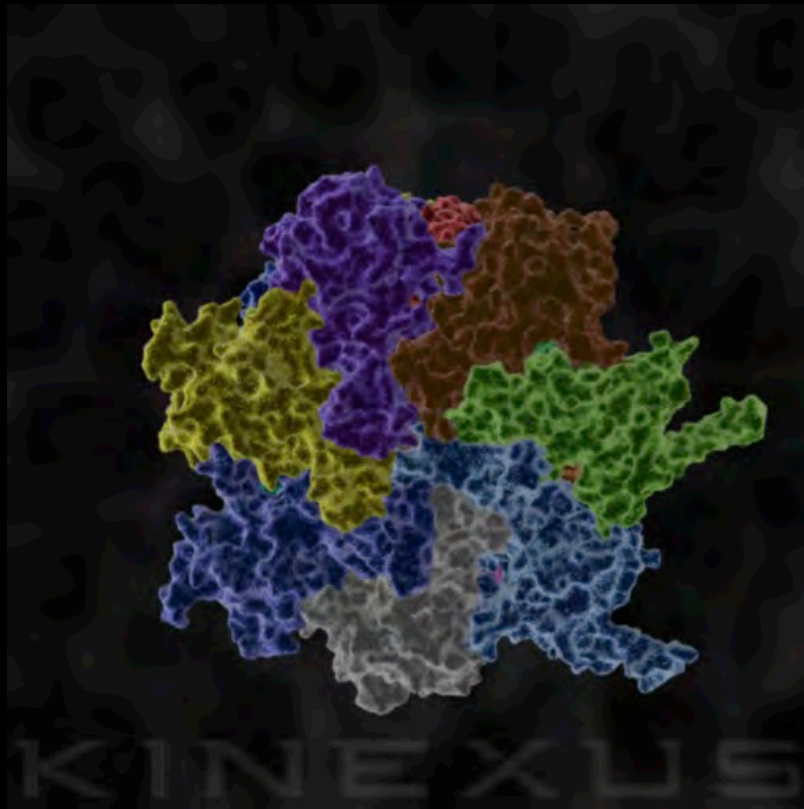


- Hyper-phosphorylation on serine and threonine may mediate the dissociation of protein complexes and protein degradation.
- Conversion to highly charged amino acids may drive protein unfolding
  - Globular proteins are converted to rods when they are coated with detergents such as sodium dodecylsulphate (SDS)
  - This is achieved because of the high negative charge on SDS; the hydrophobic part of the detergent binds to the protein and the negative charge of the sulphate permits solubilization
  - Similarly, the incorporation of highly charged clusters of phosphosites in proteins may also facilitate disruptions of their 3D structures
  - Unfolded proteins should be more susceptible to proteases

## 20S Proteasome

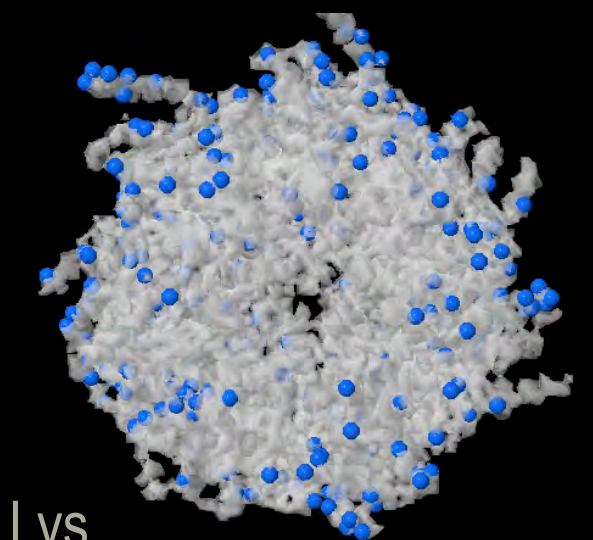
Hyperphosphorylation may facilitate protein degradation

- 20S proteasome mediates protein degradation into amino acids

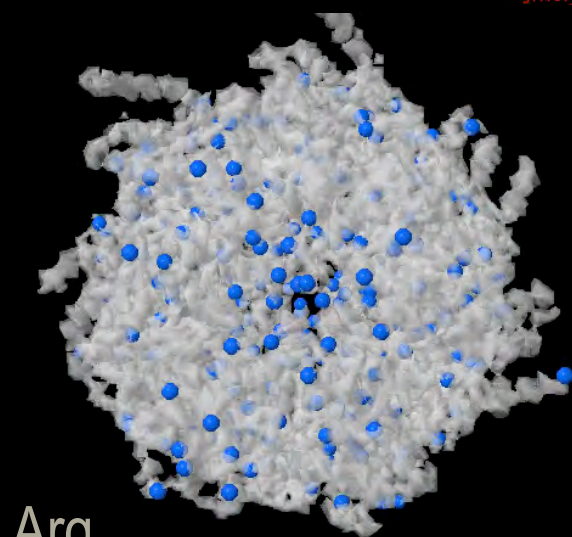




## 20S Proteasome Active Site

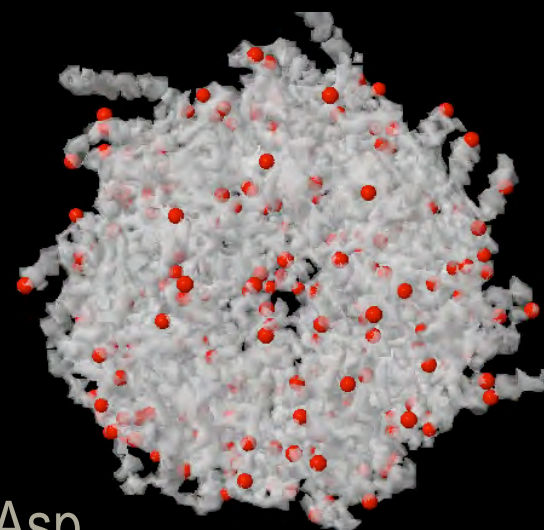


Jmol\_S

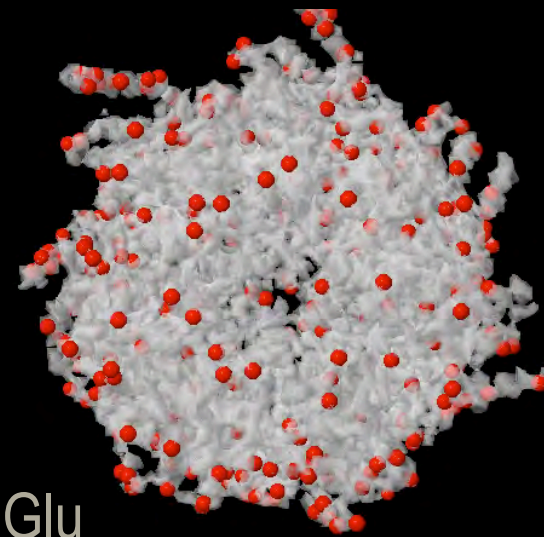


Jmol\_S

Short Chain



Jmol\_S



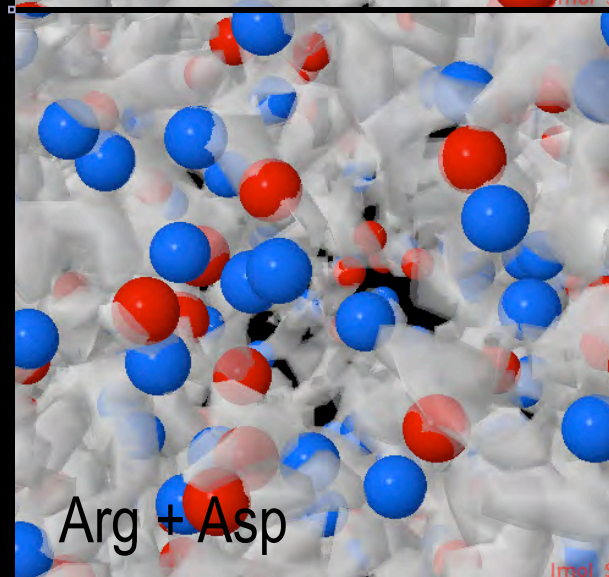
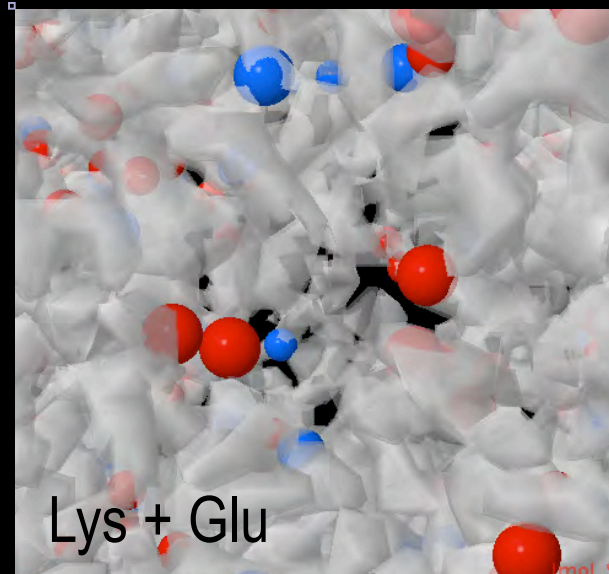
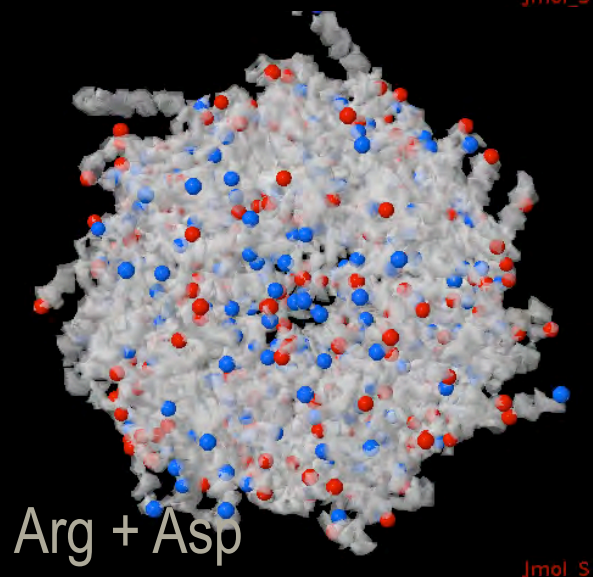
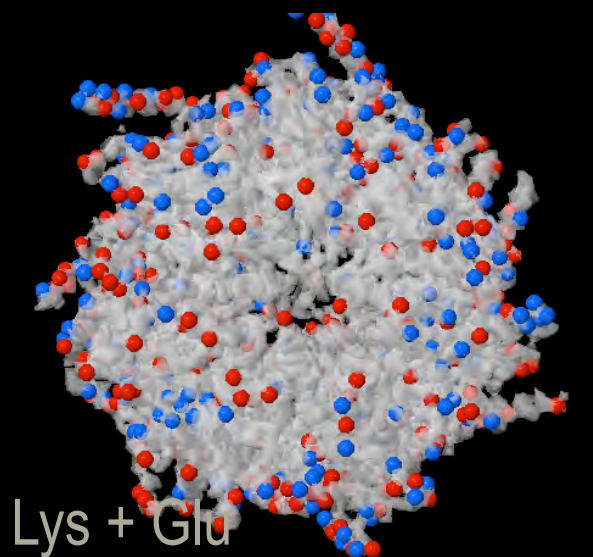
Jmol\_S

Long Chain

Glu

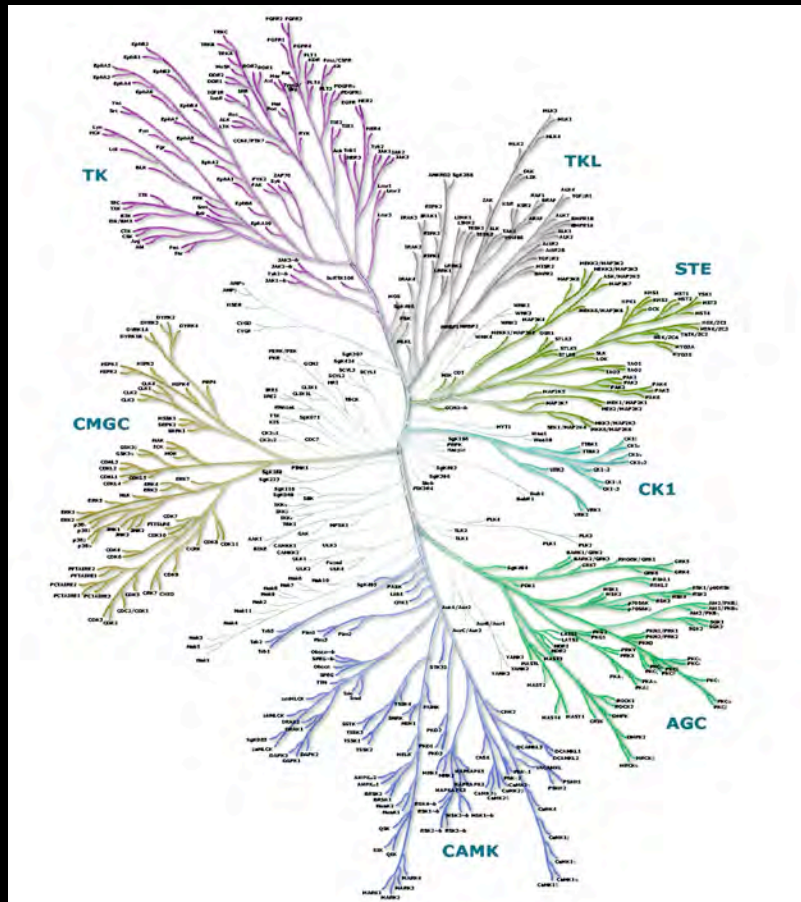


## 20S Proteasome Active Site





### The Source of the First Protein Kinase Gene



From Manning *et al.* Science 6 December 2002

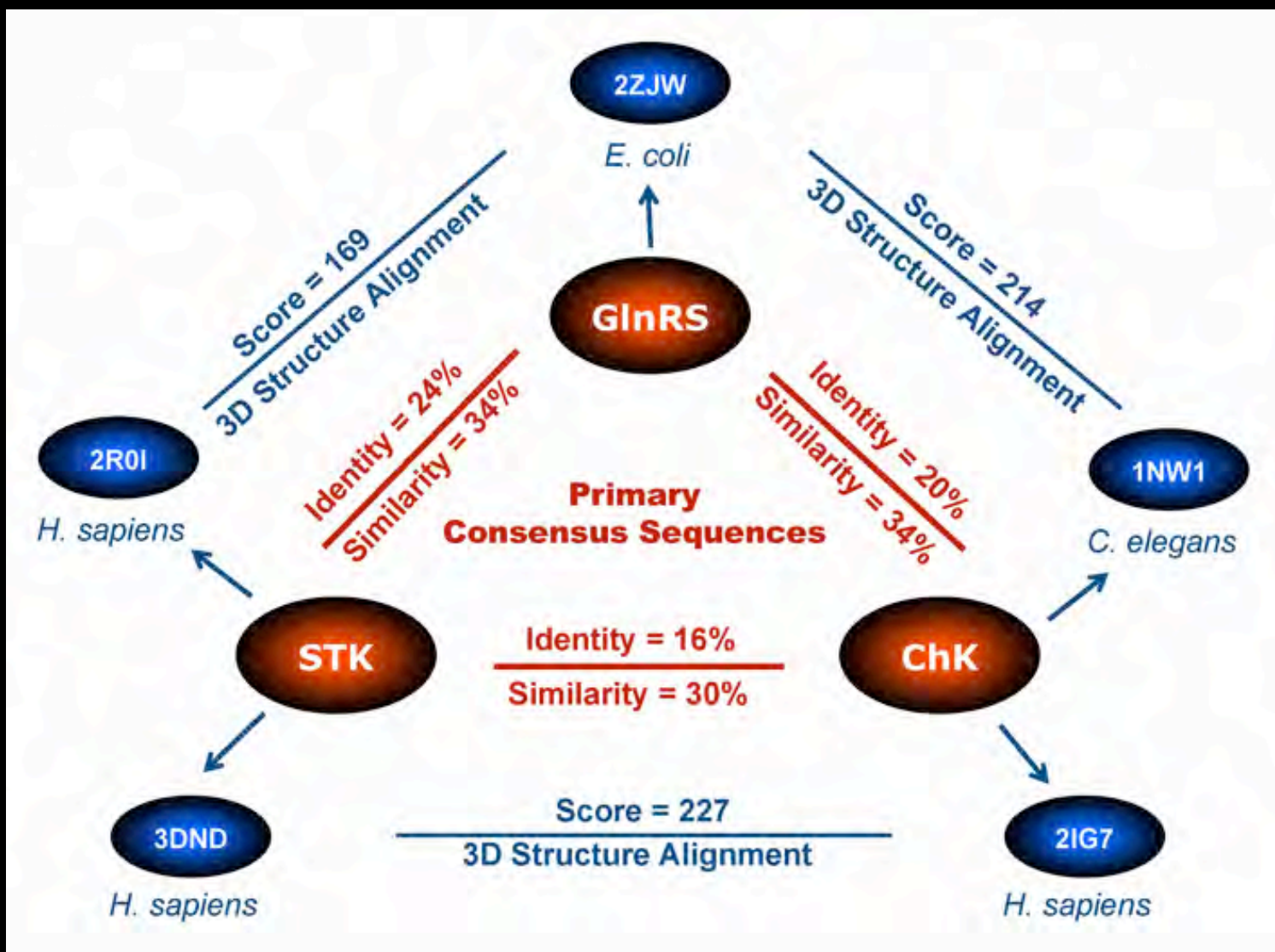
- Aligned catalytic domains of protein-serine/threonine kinases to derived consensus sequence with 247 aa
- Blast other species and ignore protein kinase sequences
- Got matches with glutamine tRNA ligase (GlnRS; for protein synthesis) and choline kinase (ChK; for phosphatidylcholine biosynthesis)
- GlnRS appears to be the source of the primordial protein kinase



# Alignment of Protein Consensus Sequences

- GlnRS = Glutamine tRNA synthetase consensus sequence
- STK = Protein-serine/threonine kinase consensus sequence
- ChK = Choline kinase consensus sequence

## Alignment of Protein Consensus Sequences



## The Emergence of Eukaryotic Cells

- Protein kinases are critical for the development of Molecular Intelligence systems in eukaryotic cells - with larger cells, communication becomes increasingly critical
- Choline kinase catalyzes the phosphorylation of choline, which is the first step in the synthesis of phosphatidylcholine and the main phospholipid found in cell eukaryotic cell membranes
- The duplication of mutation of the glutamine tRNA ligase gene on two separate occasions may have been major steps for the development of successful eukaryotic cells



- Shenshen Lai  
Ph.D. Student  
UBC



- Dr. Dirk Winkler  
Lab Manager,  
Kinexus



- Javad Safaei  
Ph.D. Student  
UBC

- Dr. Steven Pelech  
President & CSO  
Phone: 604-323-2547 Ext. 10  
E-mail: [spelech@kinexus.ca](mailto:spelech@kinexus.ca)

