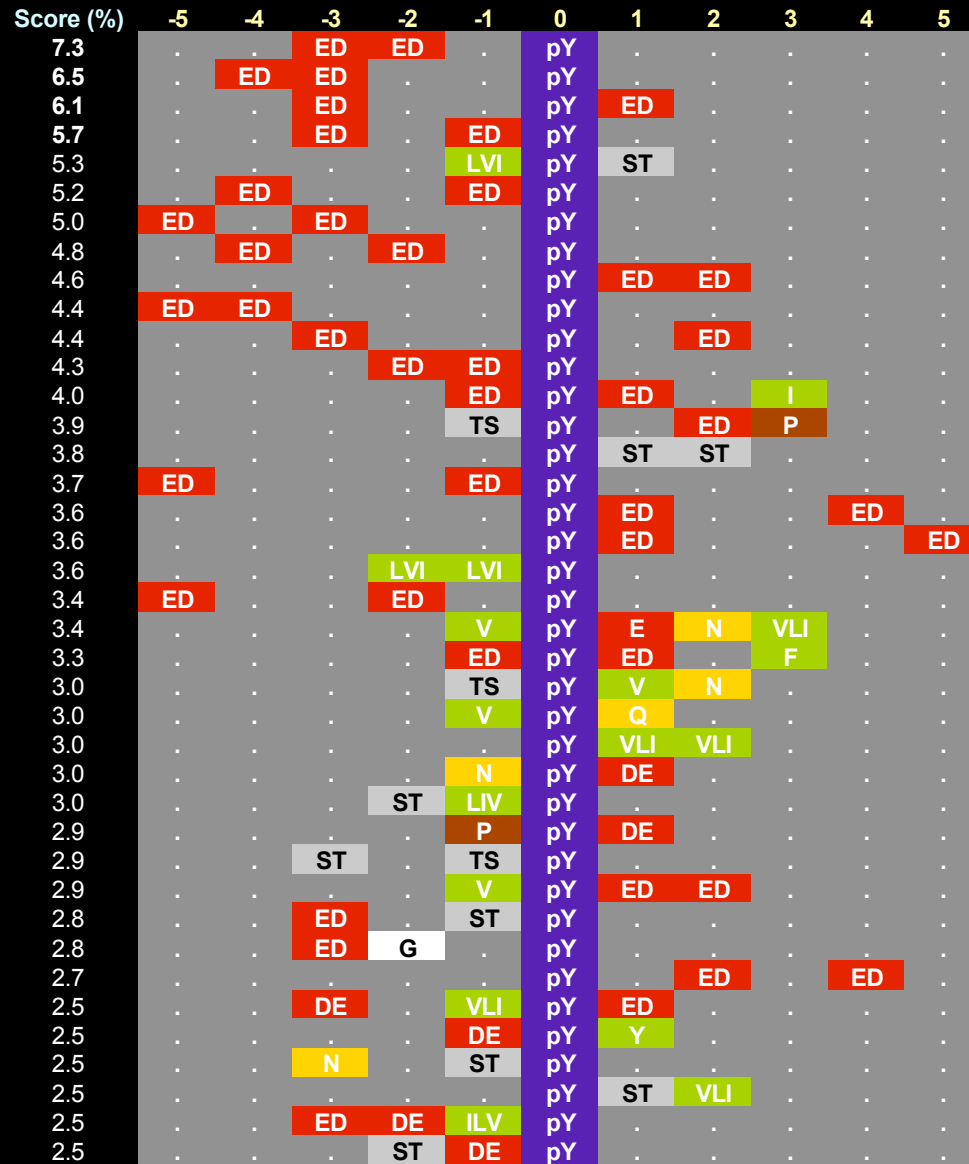


# Human Phosphoproteome - Pho-Tyr Site Frequency



Acidic	DE	Asp/Glu
Basic	KR	Lys/Arg
Weak Basic	NQ	Asn/Gln
Weak polar	M	Met
Hydrophobic	FILVY	Phe/Ile/Leu/Val/Tyr
Kink	P	Pro
Polar	ST	Ser/Thr
Phospho-AA	pSpT	pSer/pThr
Phospho-AA	pY	pTyr
Neutral	AG	Ala/Gly
Any AA	.	Variable

