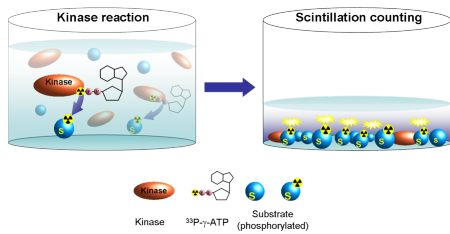


Introduction

Gene expression is regulated by multiple mechanisms. The structure of the chromatin has been identified as one major regulator of gene expression. Histone proteins constitute the core of the nucleosomal particles, which are essential for structuring the DNA into chromatin. Various post-translational modifications of histones mainly in the N-terminal region of these proteins are compacting or opening the chromatin which in turn changes the accessibility of the DNA. Such histone modifications like methylation, acetylation and phosphorylation can therefore result either in activation or in repression of the transcription of the respective genes, depending on the specific amino acid residues modified. While histone-acetylation and -methylation are currently in the focus of epigenetic research, it is also well documented that histones are extensively modified and regulated by phosphorylation. In this study we used the ProQinase *KinaseFinder* screening approach to identify kinases able to phosphorylate the N-termini of human core histone proteins under in-vitro conditions.

Principle of the assay

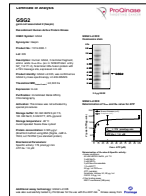


A radiometric, filter-plate based assay setup was used to determine the phosphorylation of Histone N-termini by 339 protein kinases. Kinase and potential substrate were incubated in presence of ATP containing ³³P-γ-ATP as tracer. After the kinase reaction was stopped, the proteins were immobilised on the filter membrane and the incorporated radioactivity measured by scintillation counting.

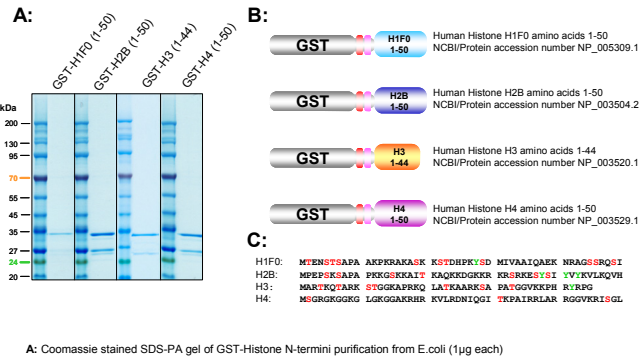
Material

Recombinant human protein kinases

- expressed in insect cells or E.coli
- tag-free or N-terminally GST- or HIS-tagged
- full length or fragments
- in-vitro activity validated for all kinases with various substrates
- details for each kinase are available at <http://proqinase.com/content/view/26>

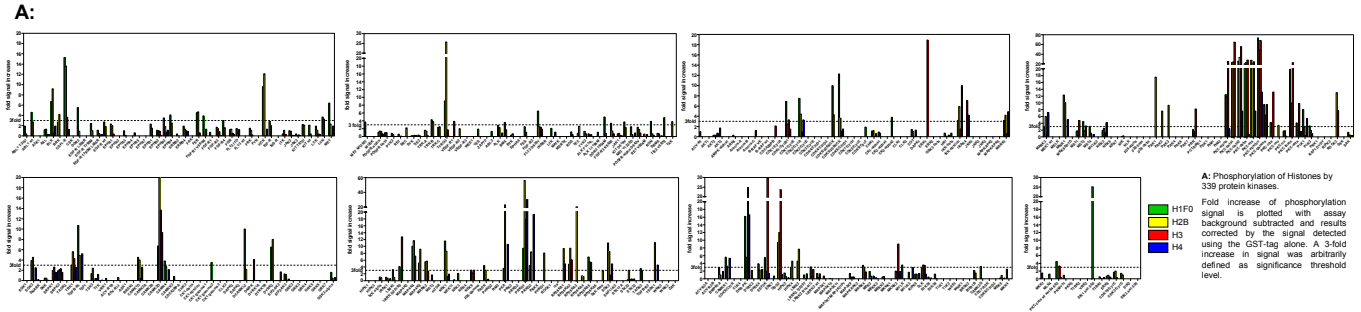


Histone substrate proteins



Results

KinaseFinder in-vitro screen of Histone phosphorylation using 339 protein kinases



B: Heatmap showing phosphorylation results for 339 protein kinases across Histone N-termini H1F0, H2B, H3, and H4. Values represent fold increase of phosphorylation signal. Green indicates significant phosphorylation (fold increase > 3).

Kinase	H1F0	H2B	H3	H4
ABL1 T319I	1.0	1.0	1.0	1.0
ABL2	1.3	0.0	0.0	0.0
ABL3	1.0	0.0	0.0	0.0
ABL4	1.0	0.0	0.0	0.0
ABL5	1.0	0.0	0.0	0.0
ABL6	1.0	0.0	0.0	0.0
ABL7	1.0	0.0	0.0	0.0
ABL8	1.0	0.0	0.0	0.0
ABL9	1.0	0.0	0.0	0.0
ABL10	1.0	0.0	0.0	0.0
ABL11	1.0	0.0	0.0	0.0
ABL12	1.0	0.0	0.0	0.0
ABL13	1.0	0.0	0.0	0.0
ABL14	1.0	0.0	0.0	0.0
ABL15	1.0	0.0	0.0	0.0
ABL16	1.0	0.0	0.0	0.0
ABL17	1.0	0.0	0.0	0.0
ABL18	1.0	0.0	0.0	0.0
ABL19	1.0	0.0	0.0	0.0
ABL20	1.0	0.0	0.0	0.0
ABL21	1.0	0.0	0.0	0.0
ABL22	1.0	0.0	0.0	0.0
ABL23	1.0	0.0	0.0	0.0
ABL24	1.0	0.0	0.0	0.0
ABL25	1.0	0.0	0.0	0.0
ABL26	1.0	0.0	0.0	0.0
ABL27	1.0	0.0	0.0	0.0
ABL28	1.0	0.0	0.0	0.0
ABL29	1.0	0.0	0.0	0.0
ABL30	1.0	0.0	0.0	0.0
ABL31	1.0	0.0	0.0	0.0
ABL32	1.0	0.0	0.0	0.0
ABL33	1.0	0.0	0.0	0.0
ABL34	1.0	0.0	0.0	0.0
ABL35	1.0	0.0	0.0	0.0
ABL36	1.0	0.0	0.0	0.0
ABL37	1.0	0.0	0.0	0.0
ABL38	1.0	0.0	0.0	0.0
ABL39	1.0	0.0	0.0	0.0
ABL40	1.0	0.0	0.0	0.0
ABL41	1.0	0.0	0.0	0.0
ABL42	1.0	0.0	0.0	0.0
ABL43	1.0	0.0	0.0	0.0
ABL44	1.0	0.0	0.0	0.0
ABL45	1.0	0.0	0.0	0.0
ABL46	1.0	0.0	0.0	0.0
ABL47	1.0	0.0	0.0	0.0
ABL48	1.0	0.0	0.0	0.0
ABL49	1.0	0.0	0.0	0.0
ABL50	1.0	0.0	0.0	0.0
ABL51	1.0	0.0	0.0	0.0
ABL52	1.0	0.0	0.0	0.0
ABL53	1.0	0.0	0.0	0.0
ABL54	1.0	0.0	0.0	0.0
ABL55	1.0	0.0	0.0	0.0
ABL56	1.0	0.0	0.0	0.0
ABL57	1.0	0.0	0.0	0.0
ABL58	1.0	0.0	0.0	0.0
ABL59	1.0	0.0	0.0	0.0
ABL60	1.0	0.0	0.0	0.0
ABL61	1.0	0.0	0.0	0.0
ABL62	1.0	0.0	0.0	0.0
ABL63	1.0	0.0	0.0	0.0
ABL64	1.0	0.0	0.0	0.0
ABL65	1.0	0.0	0.0	0.0
ABL66	1.0	0.0	0.0	0.0
ABL67	1.0	0.0	0.0	0.0
ABL68	1.0	0.0	0.0	0.0
ABL69	1.0	0.0	0.0	0.0
ABL70	1.0	0.0	0.0	0.0
ABL71	1.0	0.0	0.0	0.0
ABL72	1.0	0.0	0.0	0.0
ABL73	1.0	0.0	0.0	0.0
ABL74	1.0	0.0	0.0	0.0
ABL75	1.0	0.0	0.0	0.0
ABL76	1.0	0.0	0.0	0.0
ABL77	1.0	0.0	0.0	0.0
ABL78	1.0	0.0	0.0	0.0
ABL79	1.0	0.0	0.0	0.0
ABL80	1.0	0.0	0.0	0.0
ABL81	1.0	0.0	0.0	0.0
ABL82	1.0	0.0	0.0	0.0
ABL83	1.0	0.0	0.0	0.0
ABL84	1.0	0.0	0.0	0.0
ABL85	1.0	0.0	0.0	0.0
ABL86	1.0	0.0	0.0	0.0
ABL87	1.0	0.0	0.0	0.0
ABL88	1.0	0.0	0.0	0.0
ABL89	1.0	0.0	0.0	0.0
ABL90	1.0	0.0	0.0	0.0
ABL91	1.0	0.0	0.0	0.0
ABL92	1.0	0.0	0.0	0.0
ABL93	1.0	0.0	0.0	0.0
ABL94	1.0	0.0	0.0	0.0
ABL95	1.0	0.0	0.0	0.0
ABL96	1.0	0.0	0.0	0.0
ABL97	1.0	0.0	0.0	0.0
ABL98	1.0	0.0	0.0	0.0
ABL99	1.0	0.0	0.0	0.0
ABL100	1.0	0.0	0.0	0.0

Summary

- significant phosphorylation of Histone N-termini was observed by multiple protein kinases
- of 339 protein kinases tested, 140 (41.3%) showed significant phosphorylation of at least one Histone N-terminus
- 58 kinases (17.1%) phosphorylated 1 Histone, 37 kinases (10.9%) 2 Histones, 33 kinases (9.7%) 3 Histones and 12 kinases (3.5%) phosphorylated all 4 Histones
- multiple kinases with documented¹ Histone-phosphorylation activity were confirmed in this study
- several kinase hits are of arguable physiological relevance due to different subcellular localisation of kinase and substrate (e.g. receptor tyrosine kinases)
- using the results of this primary screen may allow to elucidate the relevance of so far unknown histone kinases