

### Abstract

Signal Transducer and Activator of Transcription 3 (STAT3) mediates cellular responses to diverse cytokines and growth factors by modulating the expression of specific target genes. While phosphorylation of STAT3 at Tyr-705 has been demonstrated to be a prerequisite for STAT3 dimerization, nuclear translocation and activation of gene transcription, the role of Ser-727 in regulation of STAT3 activity is controversial. Using Kinetworks™ KPSS 1.1 phospho-site immunoblotting screening of nocodazole-treated HeLa cells. STAT3 Ser-727 phosphorylation was enhanced during mitosis, and this correlated with a reduction of Tyr-705 phosphorylation. Over-expression of STAT3 mutants in which these phosphorylation sites were separately abolished revealed that phosphorylation at these sites appeared to be mutually antagonistic. The nocodazole-induced STAT3 Ser-727 phosphorylation was reduced by selective inhibition of cyclin-dependent kinase 1 (CDK1) phosphotransferase activity, and CDK1 could directly phosphorylate GST-STAT3 Ser-727 in vitro and it coimmunoprecipitated with STAT3 from cell lysates. Blocking Ser-727 phosphorylation through transfection of HEK-293 cells with its nonphosphorylatable mutant enhanced STAT3 DNA binding activity towards its target gene promoters, implying a negative effect of Ser-727 phosphorylation on its transcriptional activity. Interference of Ser-727 phosphorylation resulted in an exit from mitotic arrest induced by nocodazole treatment and a cell cycle arrest at G1 phase, as indicated by the accumulation of 2N cell population and enhanced expression of G1 cell cycle regulators including p21 april p27 cyclin E. Our observations point to a novel role of STAT3 Ser-727 phosphorylation in control of the onset and maintenance of M-phase during the cell cycle through downregulation of CDK inhibitors

### Introduction

STAT3 belongs to a family of signal transducing transgription factors that play gritical roles in mediating cellular SIAI3 beingstio a tamiy of signal tanksiung tamisription techs that play unital roles in med ating evaluate responses to diverse of yolines and growth factors by modeling the expression of spacific target grows. phosphorylation effort by receptor-associated tynasine kinases such as in the Janue kinase (JAK) and 50 times families or by intrinsir receptor-phorose kinases such as the explorer allowable factor response HC (EGFR). This leads to STAI3 dimetation through reciprocal phosphoryloxies-BH2 domain interactions, nuclear translocation and binding te enhance/princed regulations and the gravity and gravity effects.

Constitutive activation of STAT3 is found in a wide variety of malionancies including cancers, and the mechanisms underlying the activation of STATs have been the focus of extensive studies in recert years. While the tyrosine phosphorylation of STATs mediated by nonreceptor or receptor tyrosine kinases seems to be the the tyrosine phosphorylation of STATs mediated by nonreceptor or neorptor tyrosine kinases series to be the panding in GSTAT advantion utilized by a large number of cyclicies and growing factos, there is also the phosphorylation of a series residue within the CCDM Huminial tries or plotion advice for domain or inned STATs. STAT2 and 6. and consegorates base - 275 (277) in STAT3 in this base net domains that the triangent ploting advice the protochargent ploting and the series of the protochargent ploting and the series of the series of the series of the series of the protochargent triangent is made series phosphorylation the direction (or indirect) fails and the series (Markes), protochargent ploting advice the advice ploting series of the advice ploting series of the advice ploting series of the series of the series of the series of the advice ploting series of the series of the advice ploting series of the advice the advice series of the series of the advice the advice been controversia

During experiments in which Kinetworks™ KPSS phospho-site profiling was performed on porodazide treated HeLa cells blocked in M phase, a marked enhancement of STAT3 S727 phosphorylation was observed. The objective of this study was to determine how this was achieved and the physiological role for this evert...

## Materials and Methods

Cell Lines and Culture Conditions - HeLa (human cervical cardinoma cells) and HEK293 (human embryonic kidney cells) were maintained in Dulbecco's modified Eagle medium supplemented with 10% fetal bovine serum penicillin (100 U/m) and streptomycin (100 uo/ml), in a humidified atmosphere containing 5% CO<sub>2</sub> The culture dishes used for HEK293 cells were treated with 2% poly-L4ysine for 20 min prior to each use

Plasmids - pGEX plasmids containing glutathione-S-transferase (GST)-STAT3 wild-type (WT) and STAT3 point mutations at S727 (replaced by A, S727A) or Y705 (replaced by F, Y705F) were constructed. STAT3 expression plasmids for mammalian cell transfection experiments were prepared by inserting the *EcoRI-Xhol* fragments containing amino acids 28-770 of STAT3 from GST-STAT3 WT or mutants into pcDNA3. Flag sequences were incorporated into their Ctermini as tags

Transient Transfection - At 40-60% confluency, HEK293 cells were transiently transfected with plasmids as indicated using Lipofectamine transfection reagent.

Electrophoretic mobility shift assays - Twenty up of total HEK293 cell lysates were used in each assay carried out using STAT3 gel shift oligonucleotide pre-labeled with [y2P]ATP. The resulting protein/DNA complexes were immunoprecipitated using anti-Flag antibody, and then resolved on a 13% native polyaorylamide gel prio to detection with a fluorescence imaging analyzer.

K/networks<sup>™</sup> Probornics Analysis - Total cell lysates were propared by sonication for 15 sec in lysis buffer [pH 7.4, 20 mM Tris, 20 mM Jp-glycerophosphate, 150 mM NaC), 3 mM EDTA, 3 mM EGTA, 1 mM Na<sub>2</sub>VO<sub>2</sub> D5% Normide P-40, and 1 mM dith/divisiol) supplemental with protease inhibitors [1 mM phenylmethanesulfonylfluoride (PMSF), 2 µg/ml

leupeptin, 4 µg/ml aprotinin and 1 µg/ml pepstatin A] and centrifugation at 100,000 rpm for 30 min at 4°C. Protein concentration was determined by the Bradford assay. The Kinetworks<sup>100</sup> KPSS 1.1 phospho-site and ysis was performed with 300 µg of total protein protocols available at www.kineusx.co). Row cytometry - Twenty-four hours after transfection with STAT3 WT or S727A HEK293 cell cultures were or

plated in the new finishes to prevent overconfluence. Nonodazie treatment was administrated at 24 h after-plating for 20 h. Followingtrypsinization, cells were fixed in cold methanol and stained with propidum iodide in the presence of RNsee A. The DNA fluorescence was measured using a FACScan; data acquisition and addisis was performed with the Californet Mark to make Andrés. analysis were performed with the CellQuest software from Becton Dickinsor

BrdU incorporation assays and immundluorescence microscopy — STAT3 WT and S727A-transfected HE/233 cells were grown on coversitips and treated with nocodazole for 20 h. Following removal of the nocodazole, the cells were then incubated with 20 kM BrdU of 6 h. Following factorian or permeabilization, the cells were stained with anti-BrdU and anti-Flao antibodies for 2 h at room temperature. Their respective signals were detected by Alexa488-labeled anti-mouse IgG and Alexa569-labeled anti-rabbit IgG secondary antibodies for 1 h in the dark. Nuclei were counter-stained with propid um iodide following RNaseA digestion.

#### Results

STATS S727 phosphorylation is an mitosis-associated event. Exposure of proliferation Hel a cells to 200 ngml nocodazole for 20 h induced a cell-cycle arrest in mtosis and this was accompanied by changes of a wide range of cell signaling proteins in their phosphorylation states, as revealed by the Kinetworks™ KPSS 1.1 phosphoprotein screen (Fig. 1) This screen utilizes phospho-site-specific antibodies to systematically track prosphored access (Fig. ). In screen unices prosphored peaks announce to systemically tasks changes in prosphorylation states of 30 known signaling proteins. Unexpected y, the phosphorylation of 5727 in STAT3 (Band 35) was enhanced by about 3.5-fold after 20 h incubation with nocodazole. Treatment of Hel a with other microtubule-interfering agents such as taxid vinblastine and colchicine at concentrations effective in casing the miticabate-mining agends address taxa, ministrane, and contracte at contract and contract and the effective in casing the mitica areast also led to the increase of STAT3 S727 phosphorylation to different degrees (Fig2A). Furthermore, increased phosphorylation of STAT3 S727 phosphorylation occurred during normal mitosis in HeLa cells synchronized by releasing serum-starved cells from Go/G1 block with serun addition (Fig. 2C).

KPSS 1.1 Phospho-sites tracked:

1. Adducino S724 2. Adducino S693 3. B23 - S4 4. CDN1 - Y15 5. CREB - S133 6. ERX1 - T202+7204 7. ERX2 - T185+Y187 9. GSK3a - Y279 9. GSK3a - X279 9. GSK3a - X279 11. GSK3B - X21 12. SAMA - Y216 11. GSK3B - Y216 11. GSK3B - S17+S221 13. John - S71 14. MEX1 - S271+S221 15. MEX16 - S1895207 16. MSK1 - S376 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 18. JSB6 19. JSB6 19. JSB6 19. JSB6 19. JSB6 19. JSB7 19. JS	19. PKBa-T308 20. PKBa-K73 21. PKCa-K534. 22. PKCa-S57 23. PKCa-S57 24. PKCa-S719 25. PKR-T451 25. PKR-T451 26. Raft-S789 27. Raft-S789 28. Raft-S789 28. Raft-S789 28. Raft-S789 28. Raft-S789 29. Raft-S789 29. Raft-S789 29. Raft-S789 29. Raft-S789 20. Ra	$\begin{array}{c c c c c c c c c c c c c c c c c c c $
--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	--------------------------------------------------------

Figure 1. Kinetworks\*\* KPSS-1.1 phospho-site profile of HeLa cells in response to 200 ng/ml azole treatment for 20 h.





Figure 2. Phosphorylation of STAT3 S727 is associated with mitosis.

Reciprocal inhibition between S727 and Y705 phosphorylation - Concomitant to the increase of STAT3 S727 phosphorylation with treatment of HeLa cells with moodaacel for 20 h, there was a reduction of STAT3 phosphorylation at Y705 (Fig. 3). We examined thet wo phosphorylation events in HEX293 cells expressing wild-type (WT) STAT3 or non-phosphorylatable mutants, S727A or Y705F. Upon 20 h nocodazole treatment, whereas the over-expression of STAT3 S72A mutant abrogated S727 phosphorylation as expected, a trigher level of S727 hosphorylation was observed in STAT3 YTOF-serversing cells as compared to STAT3 WTtransfected cells (Fig.3D). Conversely, Y705 phosphorylation was further enhanced in S727A-expressing cells relativet o that in WT-transfected cells.

CDK1 phosphorylates STAT3 at S727 - Several protein-serine/threonine kinase inhibitors were employed along with nocotazole to examine their effects on STAT3 S275 phosphorylation; these included the Mek1/2 inhibitor PD98059, the p38 MAPK inhibitor SB203580, the JNK1/2 inhibitor SP600125, the CK2 inhibitor DRB (5.6-dichloro-1-beta-D-ribofuranosylbenzimidazole), the PI 3-kinase inhibitor LY294002, the CDK1/2 inhibitor olomoucine, the mTOR inhibitor rapamycin, and the PKC inhibitor R0318220. Among them, only olomoucine peturbed nocodazole-induced S727 phosphorylation significantly, while no effects were observed on STAT3 protein expression, indicating the possible involvement of CDK1 in the pathway leading to S727 phosphorylation (Fig. 4 left). Similar inhibitory effect on STAT3 S727 phosphorylation was also observed in Roscovitine, another potent and selective inhibitor of CDKs (Fig. 4 right).



Figure 3. Reciprocal inhibition between \$727 and \$705 phosphorylation in \$TAT3.



#### Figure 4. CDK1 mediates nocodazole-induced STAT3 S727 phosphorylation

To confirm the role of CDK1 in STAT3 S727 phosphonylation, we tested whether STAT3 could be phosphonylated at S727 by CDK1 in witro. Purified active recombinant CDK1/cyclin B1 complex from transfected insectcells phosphonylated GSTSTAT3at5727 (Fp.5).

GST-STAT3 - pS727 GST-STAT3 Ab CDK1 - pT161 Ab CON ST



To further explore the possible regulation of STAT3 S727 phosphorylation by CDK1, we next investigated whether CDK1 directly interacted with STAT3 in vivo. Reciprocal immunoprecipitations with either CDK1 or Rag artiblody were carried out in lysates of HEX293 of Itariastected with either a Rag-tagged STA13 WT or a S727A mutant construct, followed by nocodazole treatment for 20 h (Fig. 6). The strongest interaction was found between CDK1 and STAT3 S727A, consistent with the kinetic relationship between an enzyme and a nseudo substrate



Figure 6. CDK1 and STAT3 form complexes in vivo

STAT3 S727 phosphorylation by CDK1 attenuates its DNA-binding activity - We compared the DNA binding activities of STAT3 WT and STAT3 S727A over-expressed in HEK293 towards a double-stranded oligonucleotide containing the consensus binding site for STAT3 (5-GAT CCT TCT GG AAT TCC TA G ATC-3') in electrophoretic mobility shift assays. As shown in Figure 7, while both proteins were expressed at similar levels, more STAT3 S727A proteins were found in complexes with DNA in the STAT3 immunoprecipitates than STAT3 WT, indicating that STAT3 S727A possesses stronger DNA binding activity when compared to its wild type counterpart.



Figure 7. Inhibition of \$727 phosphorylation enhances \$TAT3 DNA-binding activity.

Inhibition of \$727 phosphorylation abolishes nocodazole-induced mitotic arrest - Cell cycle analysis using flow cytometry revealed that about 40% of the cells population transfected with STATS WT accumul ated in a tetraploid (4N) state upon a 20-h nocodazole treatment (Fig.8). Examination of cell morphology as well as chromosome configuation indicated that most cells were in mitblic arest induced by nocodazele as expected. In sharp contrast, transfecting STATS S727A resulted in the accumulation of around 70% of the cells in 2N DNA states while only less than 5% of the cells remained as 4%, indicative of a cell cycle arest at G1 phase (Fig.8). The degrease of 4N cells coupled with the increase of 2N cells implies that the mitotic arrest induced by locodazd e was lost and the cells underwent a mitotic exit and proceeded into and remained in the next G1 phase



### Figure 8. Inhibition of \$727 phosphorylation results in loss of mitotic arrest induced by nocodazole and a G1 arrest in the next cell cycle.

To further examine the molecular mechanism underlying the role of \$727 phosphorylation in nocodazole-induced mitotic arrest, we evaluated the effects of \$7A73 \$727A transfection on the expression of 30 cell cycle-related pretiens using Kinterworks" KCCP-10 cell cycle preterior analysis (Fig. 39). In particular, in the STATS S7214 transferded HEK233 cells, the expression levels of CDK4, CDK5, IEK2, 14-33 and p27 CDK inhibitor were elevated greater than 1.6-fdd, and the expression levels of Cdc34, CDK7, and cyclin G1 were reduced by more than 50% (Fig. 9B). The expression-related changes of the these and other cell cycle proteins (not shown) also point to the abrogation of nocodazole-induced mitotic arrest upon the inhibition of S727 phosphorylation.



Figure 9. Kinetworks™ KCCP 1.0 cell cycle screen analysis of the effects of STAT3 S727 phosphorylation in HE K-239 cells.

If the STAT3 S727A-expressing cells were arrested in G1 phase as indicated above, then they should enter Sphase (DNA synthesis) sconer than their wild-type counterparts when they are released from nocodazole treatment. Over 60% of STAT3 S727A cells exhibited the BrdU incorporation at 6 h after nocodazole removal In contrast, less than 15% of STAT3 WT cells were BrdU-positive (data not shown)

# Conclusions

- CDK1 directly phosphorylates STAT3 S727 during mitosis to inactive this transcription factor until critical mitotic events are completed.
- 2. The loss of phosphorylation of STAT3 S727 permits the activation of STAT3 function to facilitate cell division and release from drug induced M phase arrest
- Kinetworks<sup>™</sup> analyses of nocadazole-arrested cells has revealed a multitude of phosphorylation changes in known target phosphoproteins and cross-reactive phosphoproteins. We recently identified one of these proteins as nucleophosmin/B23. and demonstrated that polo-like kinase 3 phosphorylation of serine-4 of B23 was critical for centromere duplication. We plan to continue to identify additional phosphoproteins that are affected during mitosis to establish their interconnection in signaling pathways and their functional roles in orchestrating the myriad of events that either facilitate the successful completion of mitosis or the initiation of apoptosis.