

Report

Benzo[e]pyridoindoles, novel inhibitors of the aurora kinases

Thi My-Nhung Hoang,^{1,6} Bertrand Favier,² Annie Valette,³ Caroline Barette,⁴ Chi Hung Nguyen,⁵ Laurence Lafanechère,⁴ David S. Grierson,⁵ Stéfan Dimitrov¹ and Annie Molla^{1,*}

¹INSERM: U823 Institut Albert Bonniot; Université Joseph Fourier; La Tronche, France; ²GREPI/TIMC-IMAG CNRS: UMR 5525—Université Joseph Fourier; Grenoble, France; ³CNRS: 5088 Université Paul Sabatier; Toulouse, France; ⁴CEA, DSV, iRTSV, CNRS, UMR 5168; Centre de Criblage pour molécules Bio-Actives; Grenoble, France; ⁵UMR 176 CNRS-Institut Curie; Institut Curie; Orsay, France; ⁶Faculty of Biology; Hanoi University of Sciences; VNU, Vietnam

Key words: mitosis, aurora kinase, small-molecule inhibitors, chromosomal passenger complex, pyridoindoles, mitotic slippage

Aurora kinases are serine/threonine protein kinases that are involved in cancer development and are important targets for cancer therapy. By high throughput screening of a chemical library we found that benzo[e]pyridoindole derivatives inhibited Aurora kinases. The most potent compound (compound 1) was found to be an ATP competitive inhibitor, which inhibited *in vitro* Aurora kinases at the nanomolar range. It prevented, *ex vivo*, the phosphorylation of Histone H3, induced mitosis exit without chromosome segregation, known phenomena observed upon Aurora B inactivation. This compound was also shown to affect the localization of Aurora B, since in the presence of the inhibitor the enzyme was delocalized on the whole chromosomes and remained associated with the chromatin of newly formed nuclei.

In addition, compound 1 inhibited the growth of different cell lines derived from different carcinoma. Its IC₅₀ for H358 NSCLC (Non-Small Cancer Lung Cells), the most sensitive cell line, was 145 nM. Furthermore compound 1 was found to be efficient towards multicellular tumor spheroid growth. It exhibited minimal toxicity in mice while it had some potency towards aggressive NSCLC tumors. Benzo[e]pyridoindoles represent thus a potential new lead for the development of Aurora kinase inhibitors.

Introduction

Aurora kinases are a family of serine/threonine protein kinases that play a key role in mitosis progression.^{1,2} Aurora A is found to be associated with both centrosomes and microtubules and it is required for centrosome duplication, entry into mitosis, formation of bipolar spindle and mitotic checkpoint.³⁻⁵ Aurora B exhibits typical passenger protein behavior during mitosis. Initially, the kinase associates with centromeres, and as mitosis proceeds, it relocates to the central spindle and the midbody. Aurora B is essential for chromosome condensation, kinetochore functions, spindle checkpoint activation and cytokinesis completion.^{2,6-8}

Aurora A and B are overexpressed in many cancers, including primary colon and breast cancer.^{1,9} Furthermore, the human *Aurora A* gene is localized to the 20q13 amplicon, which is associated with a poor prognosis in breast cancer.⁹ Xenografts of mouse NIH-3T3 cells overexpressing Aurora A give rise to tumors in nude mice, suggesting that Aurora A behaves as an oncogene.¹⁰ Under similar conditions, overexpression of Aurora B may induce metastasis.¹¹ In the light of these observations, Aurora kinases have emerged as druggable targets for cancer therapy and thus, identification of Aurora kinase small molecule inhibitors is of particular interest.¹²⁻¹⁴

Several Aurora A and B inhibitors, including ZM477439, Hesperadin, VX-680 (MK-0457), MLN8054, PHA-739358 have been described.¹⁵⁻²⁰ Most of them have been included in clinical trials. VX-680, considered as the Aurora reference inhibitor, suppresses tumor growth *in vivo* and encouraging results were reported for three patients with refractory Chronic Myeloid Leukemia^{21,22} perhaps partly through the inhibition of the T315I mutant BCR-ABL.²³

The *in vitro* high throughput screening of the proprietary Institut Curie-CNRS small molecule library of 6560 compounds allowed the identification of benzo[e]pyridoindoles as new inhibitors of Aurora kinases. *Ex vivo* assays in HeLa cells showed that the identified best hit (compound 1, C1) inhibited the activity of Aurora B and mimicked the phenotype obtained after siRNA suppression of Aurora B expression. Experiments with several other tumor cell lines demonstrated that the C1 significantly affected tumor cell growth in either two- or three-dimensional cultures. C1 when used in the treatment of mice bearing H358 tumors suppressed tumor cell growth. Benzo[e]pyridoindoles are thus proposed as leads for the development of Aurora kinase inhibitors.

Results

High throughput screening for identification of aurora kinase inhibitors. We have developed an Aurora kinase assay for the identification of small molecule inhibitors of Aurora kinases. The assay is based on the phosphorylation of recombinant histone H3 by a purified recombinant Aurora A kinase domain under non-saturating ATP concentrations (see Materials and Methods for detail). The calculated statistical Z' factor was 0.77 making the assay suitable for high throughput screening. In the screen of the Institut Curie-CNRS proprietary library (chimiotheque-nationale.enscm.fr), which contains 6560 distinct mono- to penta-heterocyclic compounds,

*Correspondence to: Annie Molla; UJF Site Santé- La Tronche; Institut Albert Bonniot; INSERM U823 eq 4 BP170; 38 042 Grenoble, Cedex 9 France; Phone: 334 76 54 95 74; Fax: 33476549595; Tel.: 33476549474; Fax: 33476549595; Email: annie.molla@ujf-grenoble.fr

Submitted: 01/06/09; Accepted: 01/18/09

Previously published online as a *Cell Cycle* E-publication:
<http://www.landesbioscience.com/journals/cc/article/7879>

forty molecules were found to inhibit Aurora kinase by more than 50% when assayed at a concentration of 15 μM .

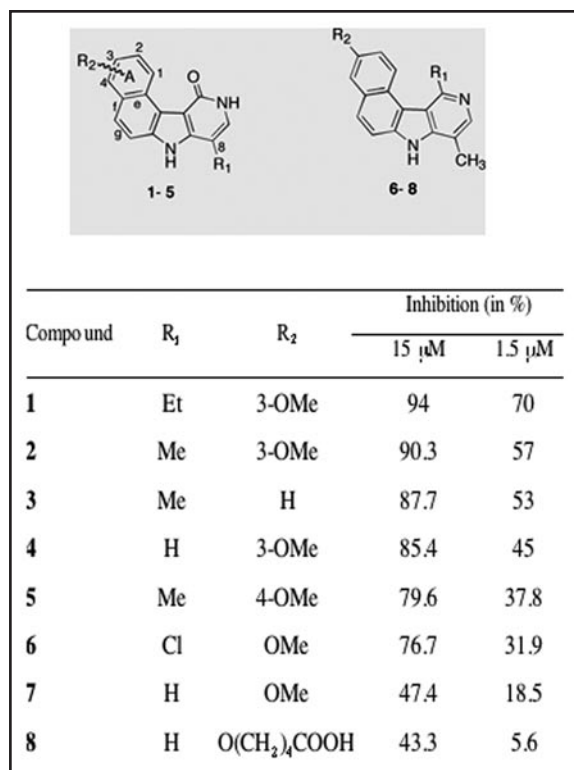
Of the fourteen most active compounds identified, six belong to the benzo[e]pyridoindole family (Fig. 1). Non-pertinent “hits” corresponding to bi- and tricyclic aromatic compounds bearing reactive (aldehyde) functionality were also found. Because of their high chemical reactivity such compounds were considered as false positive and discarded.

The most potent molecules 1 and 2 exhibit an OMe group at the 3-position of the benzo (A-ring). However, the observed high activities of compounds 3 and 5 allow the conclusion that this position of the OMe group is not a requirement for activity, and can be further moved to the 4-position (Fig. 1). The presence of an alkyl chain at position 8 is, however, crucial for the inhibitory effect of these molecules, since its absence resulted in a clear loss in activity (compare 1 with 4). Replacement of the pyridinone carbonyl oxygen by a hydrogen (7) or a chlorine atom (6) resulted also in a loss of activity. Finally, the double modification in which R1 and R2 were changed, as in 8, led to almost complete loss of activity as measured at 1.5 μM .

In an independent study (C. Cochet, personal communication) compounds 1–8 were evaluated *in vitro* for their activity against Casein Kinase II (CKII). In this screen compound 6 was found active. The other indole compounds displayed essentially no affinity for CKII. With this in mind, we have further focused on the characterization of the biological activity of benzo[e]pyridoindole 1 (C1), the compound showing the strongest inhibitory effect (Fig. 1), and we have compared its activity with that of the reference Aurora kinase inhibitor VX-680.²²

The specificity of the molecule C1 (at 1 μM concentration) towards a panel of 36 different kinases was analyzed (Fig. 2). Most of the kinases tested were not affected by the presence of C1 (Fig. 2A). C1 targets essentially Aurora A, MELK and FGFR1 kinases. These kinases were inhibited by 91, 96 and 92%, respectively (Fig. 2A). C1 had also some inhibitory activity towards CHK1, MEK1, RSK1, RSK2 but did not target the cell cycle kinases CDK and PLK (Fig. 2A). Among the Aurora kinase family, C1 exhibited an IC_{50} of 61 nM, 31 nM and 124 nM for the Aurora kinases A, B and C, respectively (Fig. 2B). These IC_{50} were similar to those determined in the same assay for VX-680, but the selectivity of the two inhibitors was slightly different. Indeed, VX-680 was more potent towards Aurora A, while the highest inhibitory effect of C1 was towards Aurora B. C1 inhibited both CHK1 (IC_{50} = 262 nM) and MELK (IC_{50} = 42 nM), whereas VX-680 was poorly efficient towards these kinases. Both inhibitors exhibited thus different selectivity (Fig. 2B), although both of them are ATP competitive Aurora inhibitors (Fig. 2C and ref. 17).

Effect of C1 on HeLa cells. The Aurora kinase inhibition efficiency of C1 was also studied *ex vivo* in HeLa cells in culture (Figs. 3 and 4). To directly demonstrate the inhibitory effect of C1 we have studied the phosphorylation status of histone H3 by western blotting (histone H3 is one of the favored substrate of Aurora B and the entry in mitosis correlates with a burst in histone H3 phosphorylation). (Figs. 3A and 4A). Treatment of the cells with C1 led to a significant decrease of H3 phosphorylation in mitotic HeLa cells (Fig. 3A). Moreover treatment of HeLa cells with compound 1 resulted in severe morphological perturbations of the HeLa cell nuclei (Fig. 3B and C).



Compound	R ₁	R ₂	Inhibition (in %)	
			15 μM	1.5 μM
1	Et	3-OMe	94	70
2	Me	3-OMe	90.3	57
3	Me	H	87.7	53
4	H	3-OMe	85.4	45
5	Me	4-OMe	79.6	37.8
6	Cl	OMe	76.7	31.9
7	H	OMe	47.4	18.5
8	H	O(CH ₂) ₄ COOH	43.3	5.6

Figure 1. Novel Aurora kinase inhibitors identified by high throughput screening. The chemical structure and the preferential conformations of the identified eight most potent hits are presented. The inhibitory efficiency (percent of inhibition) towards the kinase domain of Aurora A for each hit at 15 μM and 1.5 μM are also shown.

In particular, cells acquired enlarged lobed nuclei. Quantification showed that treatment with either C1 (2 μM) or VX-680 (300 nM) resulted in the generation of a comparable number of cells with such irregular nuclei (21% for C1 and 23% for VX-680, Fig. 3C). Note that under the same conditions, only ~1% of the control non-drug treated cells exhibited irregular nuclei (Fig. 3C).

Treatment with C1 led also to cell polyploidy (Fig. 3D). Mitotic (or 4N) cells represent about 12.5% of the population in the control, whereas they increase to 30% in the presence of compound 1 (Fig. 3D). Moreover the drug-dependent generation of enlarged nuclei (Fig. 3B) correlates with a significant increase of ploidy (58% of the cells are polyploid in the presence of 700 nM compound 1, Fig. 3D). These phenotypes (lobed nuclei and polyploid cells) were associated with an inactivation of Aurora-B,¹ suggesting that C1 is an efficient inhibitor of the Aurora B kinase activity.

Following an overnight paclitaxel treatment, the bulk mitotic cells tested positive for phospho-histone H3 (~37% of the total cell population), whereas only 3% of the total cell population was scored positive in the presence of compound 1 (2 μM) as compared to 1% for VX-680 (300 nM), the reference Aurora B kinase inhibitor (Fig. 4A and B). Note that upon C1 treatment the 4N cells number increased ~3-fold (Fig. 3D), but nonetheless a very strong inhibition of H3 phosphorylation was observed in the drug treated cells (Fig. 4A–C). The depletion of Aurora B kinase or its inactivation allowed paclitaxel treated cells to escape mitosis (mitotic slippage).¹⁶ If C1 is really responsible for inactivating Aurora B, mitotic

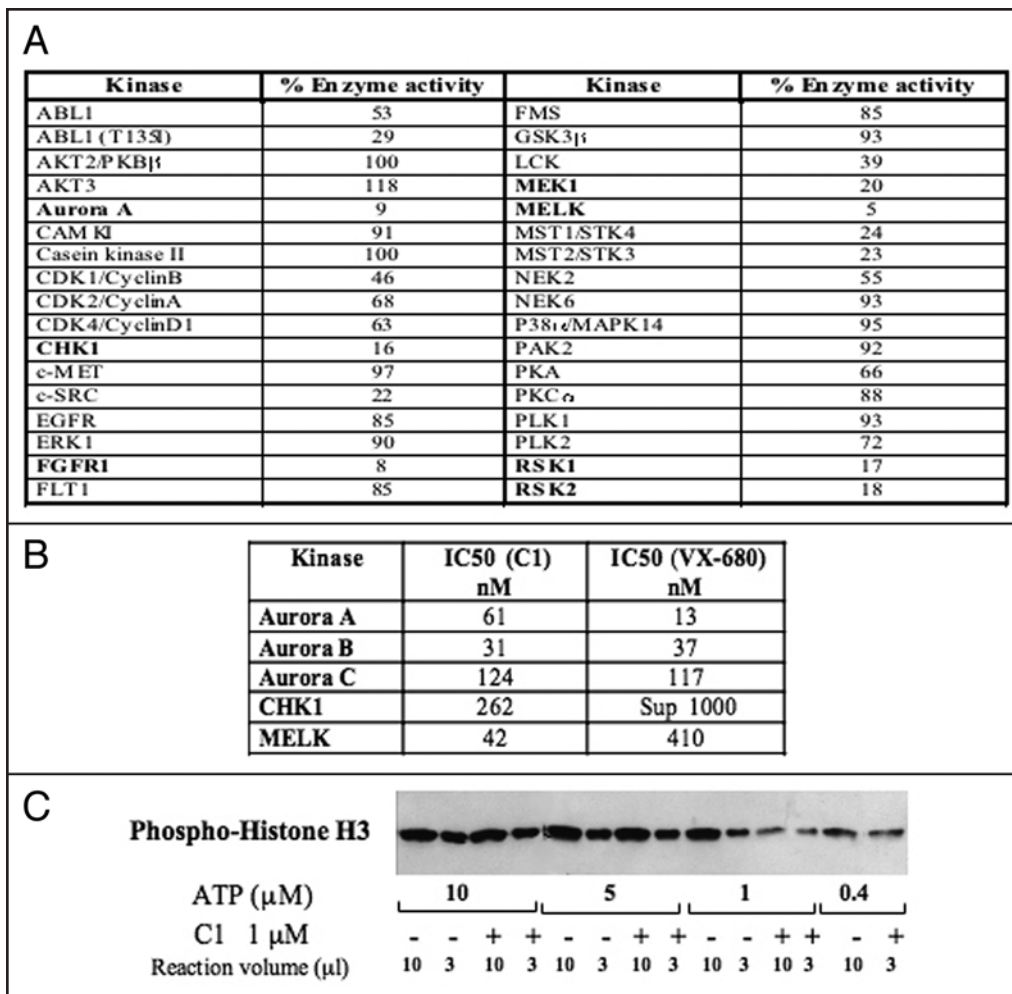


Figure 2. Kinase inhibition selectivity of C1. (A) C1 kinase profiling. 36 recombinant kinases were used in the study. The percentage of the remaining kinase activity measured upon treatment with 1 μ M C1 in a solution containing 1 μ M ATP is presented. Kinases, inhibited by more than 80%, are indicated in bold. (B) IC₅₀ values for the five best kinase targets of C1. For comparison the IC₅₀ for VX-680 for the same kinases is also shown. The measurements of IC₅₀ for both compounds were carried out under identical conditions. (C) Western blot analysis of the effect of C1 on the phosphorylation of histone H3 by Aurora A domain kinase. A kinase assay was carried out either in absence (-) or in the presence (+) of C1 (1 μ M) and at decreasing concentrations of ATP. Each reaction contained the same amount of histone H3. After completion of the reaction, 10 and 3 μ l of the reaction mixtures containing either 10, or 5 or 1 μ M ATP were run on an SDS gel (for the reaction carried out in the presence of 0.4 μ M ATP, an aliquot of only 10 μ l was run). After blotting, the phosphorylation of histone H3 was revealed by a specific antibody against the phosphorylated histone H3.

slippage should also be observed. With this in mind, we treated cells overnight with paclitaxel (33 nM), then added increasing C1 concentrations and pursued the incubation for two hours (Fig. 4C). The percentage of lobed nuclei was negligible in the control and in the 0.5 μ M C1 treated cells. The increase of C1 concentration up to 1 μ M dramatically affected the cells. Indeed, while 37% of the cells were in metaphase in the control, this percentage decreased to less than 5% in the presence of 1 μ M C1. At 0.5 μ M C1 essentially no polyploid cells and no cells with irregular nuclei were observed. At 1 μ M C1 a drastic increase of such "defective" cells was detected, i.e., ~22% of the total cell population exhibited such morphological defects (Fig. 4C). Together, these data show that cells treated with both paclitaxel and C1 exhibited mitotic slippage and therefore, reinforce the conclusion that C1 is an inhibitor of the Aurora B kinase activity.

The absence of Aurora B or the inhibition of its activity affected chromosome segregation¹⁵ and thus, treatment with C1 was expected to have the same effect, which should result in the strong polyploidy already observed (Fig. 3). To study this in real time, we have used time-lapse microscopy and stable cell lines expressing GFP-H2A, cells which allow visualization of chromatin DNA (Fig. 4D and E). The control cells (without drug treatment) showed a typical mitotic behavior with properly aligned chromosomes at metaphase, which, as mitosis proceeded, were segregated and gave finally rise to the nuclei of the two daughter cells (Fig. 4D). The treatment with 1 μ M C1 inhibited the chromosome alignment and the formation of metaphasic plate, chromosomes failed to segregate, and with time they lose their rodlike shape, decondensed and formed interphase like structure (Fig. 4E). In all analyzed cells (more than 30 in three separate experiments), the interphase type chromatin was always connected by bridges (see Fig. 4E, 46 min). These data demonstrate that upon treatment with C1 cells exit from mitosis without segregating chromosomes.

Interference with the function of the protein passenger complex resulted in a distribution of its members on the entire chromosome.² Treatment with C1 inhibited the enzymatic activity of Aurora B, and thus one should expect similar effects in the presence of C1. To test this we have followed in real time

by time-lapse microscopy the mitotic behavior of Aurora B-GFP in stable HeLa cell lines in the presence or absence of C1 (Fig. 5). Aurora B-GFP, in the control cells, had a typical passenger protein localization, decorating the centromeres at metaphase and then transferring to the central spindle and the midbody as mitosis proceeded (Fig. 5A). As expected for an Aurora B inhibitor, C1 prevented centromere alignment on the metaphasic plate (Fig. 5A and C1 500 nM). Treatment with 1 μ M C1 resulted in the partial dissociation of the GFP-tagged Aurora kinase from the centromeres (40 minutes point, Fig. 5A) and with time, Aurora B-GFP decorated also the entire chromatin (100 minutes, Fig. 5A). A similar behavior of Aurora B-GFP was observed upon treatment with VX-680, the effect being observed with a concentration of 300 nM (Fig. 5A). Immunofluorescence microscopy showed that the treatment with C1 of paclitaxel arrested mitotic cells resulted in a partial

redistribution of the endogenous Aurora B on whole chromosomes (Fig. 5B), a result in agreement with the time-lapse microscopy data.

In summary, these data illustrate in real time that the treatment of the cells with C1 affected similarly to VX-680 the behavior of Aurora B during the cell cycle, but with two-to three-fold smaller efficiency (1 μ M of C1 was necessary to observe effect close to those found with 300 nM VX-680, see Fig. 5).

Effect of C1 treatment on tumor cell line viability. The viability of tumor cell lines upon C1 treatment was investigated and compared with the effect of VX-680. Four distinct cell lines (HeLa, LL/2, H358 and HCT-116) were grown as adherent cells and the concentration that inhibits growth by two (IC_{50}) was determined (Fig. 6A). Compound 1 was found to be more potent than VX-680 to inhibit growth of HeLa cells. In the other cell lines studied it was found to be between 1.6 to 8.9 less potent than the VX-680 (Fig. 6A). Hematopoietic and lung cancer cells behaved mostly like HCT116 cells, i.e., the effect of C1 was six- to seven-fold less than this of VX-680 (data not shown). We have also investigated the effects of C1 on the HCT-116 carcinoma cells growing as MCTS by measuring the volumes of the spheroids. After treatment for 5 days the IC_{50} for HCT-116 spheroid growth inhibition was 1,066 nM for compounds 1, while it was of 108 nM for VX-680 (Fig. 6A).

Effect of C1 on tumor growth in mice. The efficiency of compound 1 was also studied in mice bearing H358 tumors. Mice bearing H358 tumors were injected with either C1 (5 mg/kg) and with the vehicle or vehicle only. The concentration of C1 was limited by its solubility in aqueous buffer. Mice were healthy after repetitive injections of the compound and grew normally. The only anomaly observed was related to the kidneys, which were slightly less colored. Blood analysis indicated neutropenia in the mice repetitively treated with C1 (average 1×10^6 neutrophil/l) compared to the control mice (average 2.6×10^6 neutrophil/l).

In the control group, tumors grew progressively and except for one, their growing index [i.e., the ratio of the tumor volume measured at the indicated time (V_D) to the initial (days 7 and 8) volume of the tumor (V_0)] was ~ 3.5 at the end of the experiment (day 51, Fig. 6B). Upon injection of C1, the tumor growth index was found to be very dispersed. Three tumors were still decreased in size at day 30, while the others behaved as control tumors. At day 37 only two tumors have regressed, while four had an index of 1 and one was growing. At day 44 two tumors increased in size (index 5.3 and 2.8) whereas two were still under control (index inferior at 1). At the end of the experiment (day 51) one tumor in the treated group has an index smaller than 1 (index 0.2). Conversely one tumor has an index superior to 6 suggesting that repetitive injections induced resistance to compound 1. Consequently the differences (mean tumor index and dispersion) between the treated and control tumors were not statistically significant (Student-t and Fisher-F test). Importantly, tumors with a decreased size (growth index below 1) were always found in the treated group after 30 days. We also

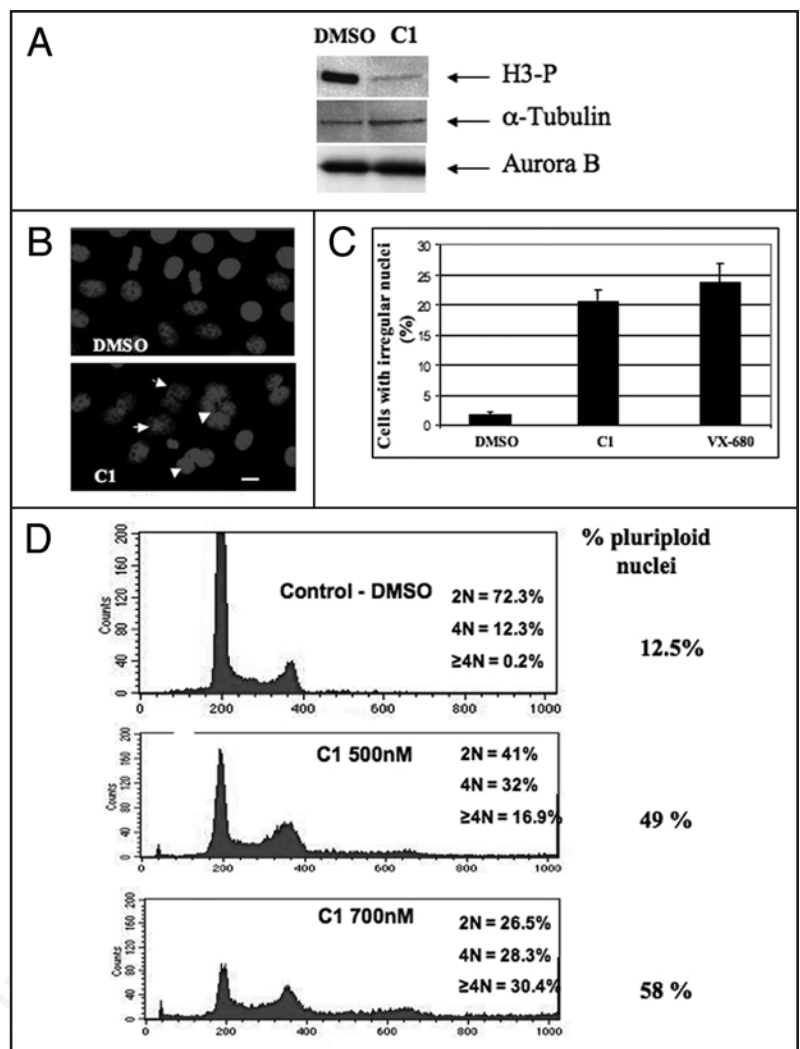


Figure 3. Effect of C1 in HeLa cells. (A) Western blot analysis of histone H3 phosphorylation. HeLa cells were incubated overnight with Nocodazole (50 nM) in the presence of either DMSO or C1 (2 μ M). Cells were then collected and lysed. Identical amounts of the lysed cell samples were run on 18% acrylamide gel containing SDS. After transfer, the blot was revealed using an antibody against phosphorylated histone H3. The same membrane was also revealed using antibodies against α -tubulin and Aurora B for estimation of both the amount of loaded proteins and of mitotic cells, respectively. (B) Treatment with C1 resulted in perturbation of the structure of the cell nuclei. HeLa cells were incubated overnight in the presence of either C1 (2 μ M) or DMSO (1%). Nuclei were stained with Hoechst 33 342 and visualized by fluorescence microscopy. Arrows indicate irregular nuclei. Similar perturbations in the nuclear structure were obtained upon treatment with 300 nM VX-680 (not shown). Bar, 5 microns. (C) Quantification of the data presented in (B). The percentage of irregular (lobed and polyploid) nuclei was determined in two independent experiments; 100 cells were analyzed per experiment. (D) FACS analysis shows that treatment with C1 results in a dramatic increase of the amount of polyploid cells. The experiments were performed with control HeLa cells (incubated only in the presence of 1% DMSO) and HeLa cells incubated with C1 (at either 500 nM or 700 nM) for 48 hours. DNA was stained with propidium iodide and the samples were analyzed by using a Beckton-Dickinson FACS analyzer. The percentage of polyploid cells (cells with a ploidy $\geq 2N$) is indicated on the right part of the figure.

found that the level of H3-phosphorylation was lower in the tumors of the C1 treated mice compared to those of the untreated control mice (Fig. 6C). Similar results were obtained in a second separate experiment (data not shown).

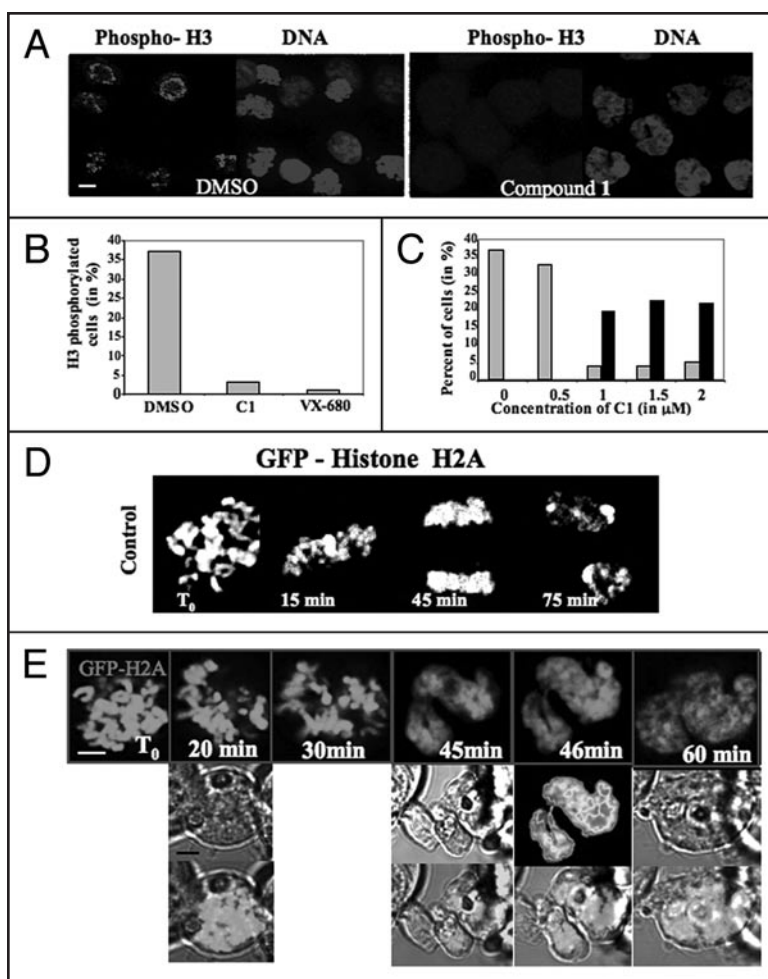


Figure 4. Effects of C1 on mitotic cells. (A) Visualization of the phosphorylation status of histone H3 in paclitaxel (33 nM) treated cells in the presence of either DMSO or C1. H3 phosphorylation was visualized by a specific antibody and DNA was stained by Hoechst 33342. Similar results were obtained with VX-680 (not shown). The bar, 5 μ m. (B) Quantification of the data presented in (A). 500 cells were analyzed in each experiment and the data of two independent experiments are shown. (C) Treatment with C1 induces mitotic slippage. HeLa cells were arrested in mitosis by overnight incubation in paclitaxel (33 nM). The cells were then incubated with the indicated concentrations of C1 for two additional hours and fixed. Mitotic and polyploid cells were scored. \sim 100 nuclei were analyzed in two different experiments; grey, mitotic cells; black, cells with polylobed nuclei. (D) Time-lapse microscopy of a stable mitotic HEK-293 cell expressing GFP-H2A histone fusion. (E) Time-lapse microscopy of C1 treated mitotic HEK-293 cells stably expressing GFP-H2A. C1 (1 μ M final concentration) was added to the overnight paclitaxel (33 nM) treated cells and then the behavior of the cell was continuously imaged. Representative photos, made at the indicated time points, are shown. The first, second and third rows present the fluorescent GFP signal and the transmission signal. The bar, 5 μ m.

Discussion

By screening a library containing 6560 mono- to pentaheterocyclic compounds, we have identified forty molecules able to significantly inhibit Aurora A kinase. The most active compounds belong to the benzo[e]pyridoindole family. These active benzo-pyridoindoles exhibit a “crescent moon” shape determined by the benzo[e] ring fusion. This geometry seems to offer the best complimentary shape to the Aurora kinase ATP site, since only two amongst the two hundred tetracyclic indole compounds screened

(including the related benzo[g] and benzo[f]pyridoindoles and linear pyrido[4,3-b]carbazoles (ellipticine analogues)) displayed some weak inhibitory activity at 1.5 μ M. The best hit (C1) was shown to be a potent in vitro, inhibitor of the three different Aurora kinases with IC_{50} of 61 nM, 31 nM and 124 nM for Aurora A, B and C, respectively. The profiling results reveal that C1 also efficiently inhibits two other kinases, FGFR and MELK. MELK, maternal embryonic leucine zipper kinase, exhibits maximal activity during mitosis but its physiological substrates are unknown.²⁴

Several different families of Aurora kinase inhibitors have now been identified and the structural basis for their inhibition selectivity was described.²⁵ For example, the tight association of either VX-680 or hesperadin with the hydrophobic pocket in the active site determined the high selectivity of the inhibitors.²⁶⁻²⁸ Note that C1, as hesperadin (a member of the SU-family of indolinone type kinase inhibitors), exhibits a cyclic amide functionality.¹⁵ Two-dimensional superimposition of C1 and hesperadin via their amide functionalities showed that the three phenyl rings in hesperadin form a “disconnected” crescent moon shape type structure, suggesting that C1 interacts with Aurora kinases in a manner similar to that of hesperadin, i.e., it may form a tight complex with the hydrophobic pocket of the kinase active site.

C1 prevents histone H3 phosphorylation in mitotic cells. This strongly indicates that in vivo the main target of C1 would be Aurora B. In agreement with this, the C1 treated cells did not exhibit monopolar phenotype, which is typically observed upon efficient inhibition of Aurora A. In addition, C1 impaired chromosome alignment and segregation as well as anaphase onset. It increased significantly the ploidy of the cells and induced mitotic slippage. All these effects were observed upon ablation of Aurora B,^{2,6} further suggesting that ex vivo the main target of C1 is this kinase. The two other strong inhibitors of Aurora kinases, VX-680 and hesperadin, were also reported to selectively target Aurora B ex vivo and to override the spindle checkpoint.^{15,28} We have also shown that both C1 and VX-680 were able to partly delocalize Aurora B kinase all over the chromosomes.

Both C1 and VX-680 efficiently prevented tumor cell proliferation. Although C1 and VX-680 have comparable in vitro inhibition efficiency towards Aurora kinases, the C1 antiproliferative activity was found to be between 1.6 and 8.9-folds weaker in the different cell lines studied compared to that of VX-680. The multicellular tumor spheroid (MCTS) model represents an intermediate level of complexity between cell growing as in vitro monolayers and solid tumors in animals. At a micromolar concentration, C1 decreased by two the spheroid volumes of HCT116 carcinoma cells growing as MCTS. The mild difference of IC_{50} in monolayer cultured cells compared to the spheroid model suggests a good biodisponibility of these compounds and its absence of multicellular resistance.

Furthermore, the effect of C1 was maximal (and comparable to that of VX-680) in H358 cell line, a cell line established from a very aggressive NSCLC tumor (Non Small Cancer Lung Cell). Patients bearing this type of tumors showed very bad prognoses and only a few relatively potent therapeutic agents against this type of tumors

are available.²⁹ With this in mind, we have further studied the *in vivo* effect of C1 on mice bearing H358 tumors. Repetitive injections of C1 were well tolerated by the animals and affected tumor growth at least for one and a half month. Note that only in few cases a decrease of the initial volume of the tumors was detected by the repetitive injections with C1 during the whole time course of the two independent experiments. This suggests that tumors escaped and became resistant to the compound upon repetitive injections.

In summary, our studies in mice revealed encouraging results with noticeable effect on a resistant model of tumors, and with few toxic effects observed so far. Benzopyridoindoles might thus be viewed as new leads for the development of Aurora kinase inhibitors that would have interesting potentials in NSCLC treatment. Combination of therapies, especially those acting via different mechanisms are proposed for improving the survival of patients with NSCLC; Aurora inhibitors, and among them Benzo[e]pyridoindole derivatives, may be included in these treatments.

Materials and Methods

Recombinant proteins. Recombinant Histone H3 and kinase domain of Aurora A were expressed in *E. coli* and purified to homogeneity. VX-680 was purchased from Kava technology Inc., while paclitaxel and nocodazole were from Sigma.

Protein kinase assay. The protein kinase assay was performed in 20 mM Tris-HCl, 20 mM KCl, 20 mM MgCl₂, 0.4 mM ATP, 0.4 mM DTT, pH 7.5. Recombinant histone H3 was used as substrate. The reaction started by the addition of the recombinant enzyme. After 1 hour of incubation at 37°C the remaining ATP was monitored by addition of kinase-Glo™ (Promega, France) under the conditions suggested by the supplier. Ten minutes later the fluorescence was recorded with a Fluostar Optima (BMG Labtechnologies). Staurosporine (0.5 mM) was used as positive control.

High throughput screening. The assay was performed as described above in black 96-well plates and started with the addition of the Aurora A kinase domain. The Z-factor of the assay was estimated to be 0.77.³⁰ The primary screening was performed in triplicate at compound concentration of 15 μM (the compounds were dissolved in 0.3% DMSO) and the selected hits were tested again at a concentration of 1.5 μM. IC₅₀ was defined as the compound concentration that leads to 50% of inhibition.

Kinase profiling and IC₅₀ determination. Assays were performed on the RBC's Discovery Dot nanoliter screening platform (Reaction Biology Corp, USA), which combines the advantages of both radioisotopes and microarrays. A kinase profiling was performed with 36 recombinant kinases. Compound 1 was tested in duplicate, at the final concentration of 1 μM. Staurosporine was used as internal control. ATP concentration was 1 μM for all reactions. The purity of the molecule 1 was superior to 95% and its synthesis is described *in*.³¹

Compound 1, VX-680 and Staurosporine were tested against five selected kinases in a 10-dose serial dilution starting at 2 μM; ATP concentration was 1 μM for all reactions. Assays were run in duplicate. IC₅₀ was defined as the compound concentration that leads to 50% of inhibition.

Cell culture. HeLa, HCT-116, Hek-293 and LL/2 were grown on Dulbecco's modified Eagle's medium (Biowhittaker, Europe). H358 cells were propagated in RPMI. Media were supplemented with 10% fetal bovine serum (Biowhittaker, Europe).

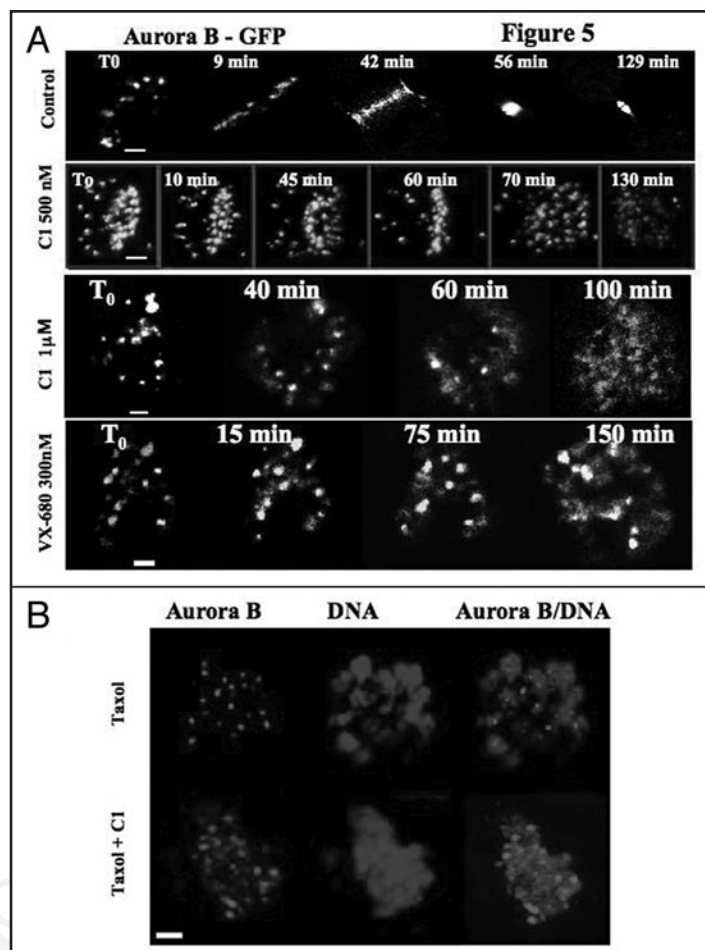


Figure 5. C1 induces the redistribution of Aurora B. (A) Time lapse microscopy of a mitotic stable HeLa cells expressing Aurora B-GFP fusion in the absence (control) or the presence of either C1 (at 500 nM or 1 μM) or VX-680 (300 nM). The compounds were added to the cell culture and then the behavior of the cells was continuously imaged. Representative photos, made at the times indicated, are presented. Note that both C1 and VX-680 delayed mitosis onset. In addition, the treatment with 1 μM C1 as well as with VX-680 resulted in a partial delocalization of Aurora B-GFP on the entire chromosomes. (B) C1 treatment induces also the redistribution of endogenous Aurora B. Overnight paclitaxel (33 nM) treated HeLa cells were incubated for two hours with C1 (1 μM), fixed and analyzed by immunofluorescence microscopy. The localization of Aurora B was detected by anti-Aurora B antibody. DNA was stained by Hoechst 33342. Note that the strictly punctuated pattern (exclusively centromeric localization) of Aurora B in the control (paclitaxel) cells is no longer observed in the C1 treated cells. Aurora B being partly diffused on chromatin.

HeLa (Aurora B-GFP) stable cell lines were already described elsewhere.³² HEK-293 cells were transfected with a plasmid expressing the fusion histone H2A—GFP and then a clonal fluorescent cell was selected and amplified.

Cell proliferation assays were conducted in 96 well cultures plates. Assays were run in triplicate. Serial dilution of compounds started at 2 μM and viability was estimated, at day 5, by addition of MTT cell counting (Promega).

The multicellular tumor spheroid (MTS) model. We have adapted the hanging-drop method³³ to produce HCT116 spheroids of similar diameters. 500 cells were suspended on the lid of an agar coated 24-petri dishes containing culture media. 48 h later the

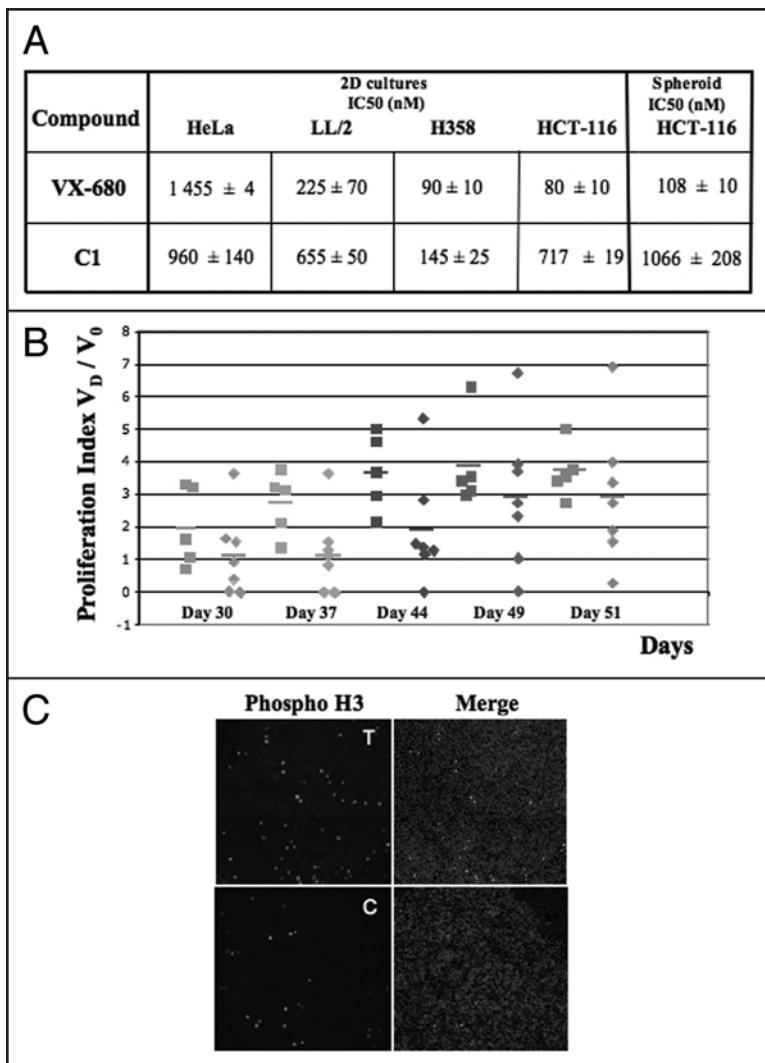


Figure 6. Effects of C1 on cell viability in two-dimensional and three-dimensional culture conditions; Effect on mice bearing H358 tumors. (A) C1 and VX-680 IC₅₀ for HeLa, LL/2, H358 and HCT-116 cells in culture (two-dimensional conditions) and for HCT-116 spheroids (3D conditions). Cell growth and viability were tested under standard conditions in 96 well culture plates with MTT (Promega) cell assay. The average of three independent experiments is shown. Note that similar IC₅₀ were determined for C1 and VX-680 towards H358 cells. (B) Proliferation indexes of H358 tumor growth. Mice were injected with 2×10^6 cells. Tumor dimensions were measured at days 7, 8, 30, 37, 44, 49 post-injection and the proliferation index (the ratio V_d/V_0 , where V_d is the volume of the tumor at the respective day, and V_0 is the average volume of the tumor calculated at day 7 and 8). Square: control mice; diamond: treated mice. Each point presents the proliferation index of one mouse tumor and the horizontal bars represent the average index of the series. (C) Histone H3 phosphorylation is decreased in C1 treated tumors. At day 51 post-injection the control and the C1 treated H358 tumor bearing cell mice were sacrificed and a small part of the tumors were immediately frozen. The phosphorylation status of histone H3 in frozen sections of the tumors of the control (C) or C1 treated mice (T) was visualized by using a specific anti-phosphorylated histone H3 antibody. DNA was stained by Hoechst 33342.

spheroids were transferred to the culture medium. Spheroid volumes were measured before drug treatment (D_0). HCT116 spheroids were treated for 5 days with different concentrations of the compounds. Control HCT116 spheroids were grown under the same conditions, but without drug treatment. The size of each spheroid was determined by measuring 2 orthogonal diameters (d_1 and d_2) using

an inverted microscope. The volume was calculated according to the formula: $V = 4/3\pi r^3$ where $r = 1/2 \sqrt{d_1 \times d_2}$. Spheroid growth was calculated by measuring the variations in volume.

Ex vivo microscopy. Ex vivo experiments were conducted on cells grown on Lab-Tek chambered coverglass (Nalge Nunc International) and maintained under standard culture conditions as described in.³² Images were acquired on a Zeiss LSM510 system using a Planapochromat 40x water immersion objective. GFP was excited with a 488 nm Argon 2 laser (power varying from 0.1 to 2%). Confocal slices are shown.

Immunofluorescence. Cells grown for 24 hours on glass coverslips were fixed at 37°C in 4% formaldehyde, 2% sucrose and then immunofluorescences were then performed as described in.³⁴ Phosphorylated Histone H3 was detected by a polyclonal rabbit antibody (Upstate). Aurora B was detected by using mouse monoclonal AIM-1 (1/100, Transduction Laboratories). DNA was visualized with 0.1 mM Hoechst 33342 (Sigma). Images were collected with a ZEISS 510 Laser Scanning Confocal microscope with a 63x immersion oil objective. Slices of 0.5 micron are shown.

Cells seeded on glass-coverslips were arrested in mitosis by the addition of paclitaxel (33 nM). Immunohistochemistry was performed on 8 nm thick frozen sections as describe above.

In vivo experiments. In vivo experiments were conducted on four-weeks old female Swiss nude mice (Iffa Credo, Marcy l'Etoile, France). After one week of adaptation in the animal facility (French agreement number A38-516-01), the mice were inoculated subcutaneously with 2×10^6 exponentially growing H358 cells mixed with growth factor free matrigel (1/1, BD). Tumors were established at seven days post-injection. Then the mice from each cage were randomly divided into two groups, which allowed the equalization of the mean tumor size of each group. Tumor volumes were determined by measuring two perpendicular diameters using a clipper and then calculated as follows: $V = d_1^2 \times d_2/2$, where d_1 and d_2 are the smallest and the largest diameters, respectively. One mice group (7 animals) received the treatment (compound 1, 100 µg per 20 g mouse in vehicle buffer (PEG 300/PBS/DMSO; 5/4/1) intraperitoneally, whereas the control group (6 mice) was injected with vehicle only. Mice were injected four times a week (Monday, Tuesday, Thursday and Friday). Once a week, mice were weighed and the volumes of the tumors were measured. Only one animal (from the control group) displayed a weight loss and was sacrificed. All other mice received 23 injections, then sacrificed and autopsied. For further analysis, blood was sampled and tumors were dissected and weighed. A piece of tumor was also embedded in OCT compound and frozen for immuno-histochemistry experiments. A second similar separate experiment that included five animals for each group was also carried out.

Acknowledgements

The authors thank Fabienne Sirot for help in the initial phase of this work and Céline Frangia for her skilful technical assistance. We thank Dr. Dimitrios Shkoufias for reading the manuscript and suggestions. Thanks to the ChemAxon company (www.chemaxon.com), which has allowed the free use of the MarvinView package. T.H. was supported by a Vietnam/French program and was a recip-

ient of an ARC fellowship. This work was supported by INSERM, ANR (Project R05075CC), ARECA and Rhône-Alpes Canceropole CLARA (EpiProNetwork). S.D. acknowledges La Ligue Nationale contre le Cancer (Equipe labellisée La Ligue).

References

- Vader G, Lens SM. The aurora kinase family in cell division and cancer. *Biochim biophys Acta* 2008; 1786:60-72.
- Ruchaud S, Carmena M, Earnshaw WC. Chromosomal passengers conducting cell division. *Nat Rev Mol Cell Biol* 2007; 8:798-812.
- Sardon T, Peset I, Petrova B, Vernos I. Dissecting the role of Aurora A during spindle assembly. *EMBO J* 2008; 27:2567-79.
- Hannak E, Kirkham M, Hyman AA, Oegema K. Aurora-A kinase is required for centrosome maturation in *Caenorhabditis elegans*. *J Cell Biol* 2001; 155:1109-16.
- Seki A, Coppinger JA, Jang CY, Yates JR, Fang G. Bora and the kinase Aurora A cooperatively activate the kinase Plk1 and control mitotic entry. *Science* 2008; 320:1655-8.
- Lens SM, Medema RH. The survivin/Aurora B complex: its role in coordinating tension and attachment. *Cell Cycle* 2003; 2:507-10.
- Ditchfield C, Johnson VL, Tighe A, et al. Aurora B couples chromosome alignment with anaphase by targeting BubR1, Mad2 and Cenp-E to kinetochores. *J Cell Biol* 2003; 161:267-80.
- Ditchfield C, Keen N, Taylor SS. The Ipl1/Aurora kinase family: methods of inhibition and functional analysis in mammalian cells. *Methods Mol Biol* 2005; 296:371-81.
- Sen S, Zhou H, White RA. A putative serine/threonine kinase encoding gene BTAK on chromosome 20q13 is amplified and overexpressed in human breast cancer cell lines. *Oncogene* 1997; 14:2195-200.
- Bischoff JR, Anderson L, Zhu Y, et al. A homologue of *Drosophila* aurora kinase is oncogenic and amplified in human colorectal cancers. *EMBO J* 1998; 17:3052-65.
- Giet R, Petretti C, Prigent C. Aurora kinases, aneuploidy and cancer, a coincidence or a real link? *Trends Cell Biol* 2005; 15:241-50.
- Keen N, Taylor S. Aurora-kinase inhibitors as anticancer agents. *Nat Rev Cancer* 2004; 4:927-36.
- Girdler F, Gascoigne KE, Evers PA, et al. Validating Aurora B as an anti-cancer drug target. *J Cell Sci* 2006; 119:3664-75.
- Jackson JR, Patrick DR, Dar MM, Huang PS. Targeted anti-mitotic therapies: can we improve on tubulin agents? *Nat Rev Cancer* 2007; 7:107-17.
- Hauf S, Cole RW, LaTerra S, et al. The small molecule Hesperadin reveals a role for Aurora B in correcting kinetochore-microtubule attachment and in maintaining the spindle assembly checkpoint. *J Cell Biol* 2003; 161:281-94.
- Gadea BB, Ruderman JV. Aurora kinase inhibitor ZM447439 blocks chromosome-induced spindle assembly, the completion of chromosome condensation, and the establishment of the spindle integrity checkpoint in *Xenopus* egg extracts. *Mol Biol Cell* 2005; 16:1305-18.
- Harrington EA, Bebbington D, Moore J, et al. VX-680, a potent and selective small-molecule inhibitor of the Aurora kinases, suppresses tumor growth in vivo. *Nat Med* 2004; 10:262-7.
- Fancelli D, Moll J, Varasi M, et al. 1,4,5,6-tetrahydropyrrolo[3,4-c]pyrazoles: identification of a potent aurora kinase inhibitor with a favorable antitumor kinase inhibition profile. *J Med Chem* 2006; 49:7247-51.
- Soncini C, Carpinelli P, Gianellini L, et al. PHA-680632, a novel Aurora kinase inhibitor with potent antitumoral activity. *Clin Cancer Res* 2006; 12:4080-9.
- Hoar K, Chakravarty A, Rabino C, et al. MLN8054, a small-molecule inhibitor of Aurora A, causes spindle pole and chromosome congression defects leading to aneuploidy. *Mol Cell Biol* 2007; 27:4513-25.
- Gautschi O, Heighway J, Mack PC, Purnell PR, Lara PN Jr, Gandara DR. Aurora kinases as anticancer drug targets. *Clin Cancer Res* 2008; 14:1639-48.
- Bain J, Plater L, Elliott M, et al. The selectivity of protein kinase inhibitors: a further update. *Biochem J* 2007; 408:297-315.
- Noronha G, Cao J, Chow CP, et al. Inhibitors of ABL and the ABL-T315I mutation. *Curr Top Med Chem* 2008; 8:905-21.
- Badouel C, Körner R, Frank-Vaillant M, Couturier A, Nigg EA, Tassan JP. M-phase MELK activity is regulated by MPF and MAPK. *Cell Cycle* 2006; 5:883-9.
- Cheetham GM, Charlton PA, Golec JM, Pollard JR. Structural basis for potent inhibition of the Aurora kinases and a T315I multi-drug resistant mutant form of Abl kinase by VX-680. *Cancer Lett* 2007; 251:323-9.
- Cheetham GM, Knegt RM, Coll JT, et al. Crystal structure of aurora-2, an oncogenic serine/threonine kinase. *J Biol Chem* 2002; 277:42419-22.
- Bayliss R, Sardon T, Vernos I, Conti E. Structural basis of Aurora-A activation by TPX2 at the mitotic spindle. *Mol Cell* 2003; 12:851-62.
- Sessa F, Mapelli M, Ciferri C, et al. Mechanism of Aurora B activation by INCENP and inhibition by hesperadin. *Mol Cell* 2005; 18:379-91.
- Alvarez M, Roman E, Santos ES, Racz LE. New targets for non-small-cell lung cancer therapy. *Expert Rev Anticancer Ther* 2007; 7:1423-37.
- Zhang JH, Chung TD, Oldenburg KR. A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. *J Biomol Screen* 1999; 4:67-73.
- Nguyen CH, Bisagni E, Lavelle F, Bissery MC, Huel C. Synthesis and antitumor properties of new 4-methyl-substituted- pyrido[4,3-b]indoles (gamma-carbolines). *Anticancer Drug Des* 1992; 7:219-33.
- Delacour-Larose MMA, Skoufias DA, Margolis RL, Dimitrov S. Distinct dynamics of Aurora B and Survivin during mitosis. *Cell Cycle* 2004; 3:1418-26.
- Del Duca D, Werbowetski T, Del Maestro RE. Spheroid preparation from hanging drops: characterization of a model of brain tumor invasion. *J Neurooncol* 2004; 67:295-303.
- Delacour-Larose M, Thi MN, Dimitrov S, Molla A. Role of survivin phosphorylation by aurora B in mitosis. *Cell Cycle* 2007; 6:1878-85.