

# SCREENING PROJECT FORM (SPF)

Please complete both parts of this form (administrative & scientific)

page 1/5  
(V18)

## Administrative Information

Project ID:  
(provided by ProQinase)

Please note: Testing can not be performed without signature (see page 5)!

Company:

Contact person:  
(appearing on report)

Address:

E-mail:

Phone:  
(optional)

Remarks:

## Scientific Information

(see our Guidelines For Biochemical Screening Services for details)

# of kinases:

# of compounds:

Single Concentration:

IC50:

(10 semi log dilutions)

Singlicate measurement:

Duplicate measurement:

(see guidelines)

### Compounds provided in:

Plates (liquid samples)

Volume of compound  
stocks (per well):

  $\mu$ l

Molarity of  
compound stocks:

 M

# of plates shipped:

Plate format:

- 96 well  
 384 well

Vials (liquid samples)

# of vials shipped:

Please provide information about  
molarity and volume of stock solutions  
using the Compound Submission Form.

Vials (solid samples)

# of vials shipped:

Please provide information about  
amount and molecular weights using  
the Compound Submission Form.

Assay concentration or highest final assay concentration (for IC50) of compounds  
will be 1/100 of stock concentration.  
(DMSO final assay concentration will be 1%)

Remarks:

## Kinases to be tested (please mark):

(list continues on following page)

### I. Protein Kinases

- |                                                   |                                           |                                               |                                           |
|---------------------------------------------------|-------------------------------------------|-----------------------------------------------|-------------------------------------------|
| <b>A</b>                                          | <b>C</b>                                  | <input type="checkbox"/> CK1-gamma2           | <input type="checkbox"/> EPHA5            |
| <input type="checkbox"/> ABL1 wt                  | <input type="checkbox"/> CAMK1D           | <input type="checkbox"/> CK1-gamma3           | <input type="checkbox"/> EPHA6            |
| <input type="checkbox"/> ABL1 G250E               | <input type="checkbox"/> CAMK2A           | <input type="checkbox"/> CK2-alpha1           | <input type="checkbox"/> EPHA7            |
| <input type="checkbox"/> ABL1 Q252H               | <input type="checkbox"/> CAMK2B           | <input type="checkbox"/> CK2-alpha2           | <input type="checkbox"/> EPHA8            |
| <input type="checkbox"/> ABL1 Y253F               | <input type="checkbox"/> CAMK2D           | <input type="checkbox"/> CLK1                 | <input type="checkbox"/> EPHB1            |
| <input type="checkbox"/> ABL1 E255K               | <input type="checkbox"/> CAMK2G           | <input type="checkbox"/> CLK2                 | <input type="checkbox"/> EPHB2            |
| <input type="checkbox"/> ABL1 T315I               | <input type="checkbox"/> CAMK4            | <input type="checkbox"/> CLK3                 | <input type="checkbox"/> EPHB3            |
| <input type="checkbox"/> ABL1 F317I               | <input type="checkbox"/> CAMK4            | <input type="checkbox"/> CLK4                 | <input type="checkbox"/> EPHB4            |
| <input type="checkbox"/> ABL1 M351I               | <input type="checkbox"/> CAMKK1           | <input type="checkbox"/> COT                  | <input type="checkbox"/> ERBB2 wt         |
| <input type="checkbox"/> ABL1 H396P               | <input type="checkbox"/> CAMKK2           | <input type="checkbox"/> CSF1-R               | <input type="checkbox"/> ERBB2 775YVMA776 |
| <input type="checkbox"/> ABL2                     | <input type="checkbox"/> CDC6-RET         | <input type="checkbox"/> CSK                  | <input type="checkbox"/> ERBB4            |
| <input type="checkbox"/> ACK1                     | <input type="checkbox"/> CDC42BPA         | <b>D</b>                                      | <input type="checkbox"/> ERK1             |
| <input type="checkbox"/> ACV-R1                   | <input type="checkbox"/> CDC42BPB         | <input type="checkbox"/> DAPK1                | <input type="checkbox"/> ERK2             |
| <input type="checkbox"/> ACV-R1B                  | <input type="checkbox"/> CDC7/DBF4        | <input type="checkbox"/> DAPK2                | <input type="checkbox"/> ERK5             |
| <input type="checkbox"/> ACV-R2A                  | <input type="checkbox"/> CDK1/CycA2       | <input type="checkbox"/> DAPK3                | <input type="checkbox"/> ERK7             |
| <input type="checkbox"/> ACV-R2B                  | <input type="checkbox"/> CDK1/CycB1       | <input type="checkbox"/> DCAMKL2              | <b>F</b>                                  |
| <input type="checkbox"/> ACV-RL1                  | <input type="checkbox"/> CDK1/CycE1       | <input type="checkbox"/> DDR2 wt              | <input type="checkbox"/> FAK aa2-1052     |
| <input type="checkbox"/> AKT1 aa1-480             | <input type="checkbox"/> CDK2/CycA2       | <input type="checkbox"/> DDR2 N456S           | <input type="checkbox"/> FER              |
| <input type="checkbox"/> AKT1 aa106-480           | <input type="checkbox"/> CDK2/CycD1       | <input type="checkbox"/> DDR2 T654M           | <input type="checkbox"/> FES              |
| <input type="checkbox"/> AKT2 aa1-481             | <input type="checkbox"/> CDK2/CycE1       | <input type="checkbox"/> DMPK                 | <input type="checkbox"/> FGF-R1 wt        |
| <input type="checkbox"/> AKT2 aa107-481           | <input type="checkbox"/> CDK3/CycC        | <input type="checkbox"/> DNA-PK               | <input type="checkbox"/> FGF-R1 V561M     |
| <input type="checkbox"/> AKT3 aa1-479             | <input type="checkbox"/> CDK3/CycE1       | <input type="checkbox"/> DYRK1A               | <input type="checkbox"/> FGF-R2           |
| <input type="checkbox"/> AKT3 aa106-479           | <input type="checkbox"/> CDK4/CycD1       | <input type="checkbox"/> DYRK1B               | <input type="checkbox"/> FGF-R3 wt        |
| <input type="checkbox"/> ALK wt (GST-HIS-tag)     | <input type="checkbox"/> CDK4/CycD2       | <input type="checkbox"/> DYRK2                | <input type="checkbox"/> FGF-R3 K650E     |
| <input type="checkbox"/> ALK C1156Y (GST-HIS-tag) | <input type="checkbox"/> CDK4/CycD3       | <input type="checkbox"/> DYRK3                | <input type="checkbox"/> FGF-R3 K650M     |
| <input type="checkbox"/> ALK F1174L (GST-HIS-tag) | <input type="checkbox"/> CDK5/p25NCK      | <input type="checkbox"/> DYRK4                | <input type="checkbox"/> FGF-R3 G697C     |
| <input type="checkbox"/> ALK F1174S (GST-HIS-tag) | <input type="checkbox"/> CDK5/p35NCK      | <b>E</b>                                      | <input type="checkbox"/> FGF-R4           |
| <input type="checkbox"/> ALK L1196M (GST-HIS-tag) | <input type="checkbox"/> CDK6/CycD1       | <input type="checkbox"/> EEF2K                | <input type="checkbox"/> FGR              |
| <input type="checkbox"/> ALK L1196M (GST-HIS-tag) | <input type="checkbox"/> CDK6/CycD2       | <input type="checkbox"/> EGF-R wt             | <input type="checkbox"/> FLT3 wt          |
| <input type="checkbox"/> ALK G1202R (GST-HIS-tag) | <input type="checkbox"/> CDK6/CycD3       | <input type="checkbox"/> EGF-R G719C          | <input type="checkbox"/> FLT3 D835Y       |
| <input type="checkbox"/> ALK R1275Q (GST-HIS-tag) | <input type="checkbox"/> CDK7/CycH/MAT1   | <input type="checkbox"/> EGF-R G719S          | <input type="checkbox"/> FLT3 ITD         |
| <input type="checkbox"/> AMPK-alpha1 aa1-550      | <input type="checkbox"/> CDK8/CycC        | <input type="checkbox"/> EGF-R G719S          | <input type="checkbox"/> FRK              |
| <input type="checkbox"/> ARK5                     | <input type="checkbox"/> CDK9/CycK        | <input type="checkbox"/> EGF-R d746-750       | <input type="checkbox"/> FYN wt           |
| <input type="checkbox"/> ASK1                     | <input type="checkbox"/> CDK9/CycT1       | <input type="checkbox"/> EGF-R d747-749 A750P | <input type="checkbox"/> FYN Y531F        |
| <input type="checkbox"/> Aurora-A                 | <input type="checkbox"/> CDK12 wt/CycK    | <input type="checkbox"/> EGF-R d747-752 P753S | <b>G</b>                                  |
| <input type="checkbox"/> Aurora-B                 | <input type="checkbox"/> CDK12 R722C/CycK | <input type="checkbox"/> EGF-R d752-759       | <input type="checkbox"/> GRK2             |
| <input type="checkbox"/> Aurora-C                 | <input type="checkbox"/> CDK13/CycK       | <input type="checkbox"/> EGF-R T790M          | <input type="checkbox"/> GRK3             |
| <input type="checkbox"/> AXL                      | <input type="checkbox"/> CDK16/CycY       | <input type="checkbox"/> EGF-R L858R          | <input type="checkbox"/> GRK4             |
| <b>B</b>                                          | <input type="checkbox"/> CDK17/p35NCK     | <input type="checkbox"/> EGF-R L861Q          | <input type="checkbox"/> GRK5             |
| <input type="checkbox"/> BLK                      | <input type="checkbox"/> CDK18/CycY       | <input type="checkbox"/> EIF2AK2              | <input type="checkbox"/> GRK6             |
| <input type="checkbox"/> Bmpr1A                   | <input type="checkbox"/> CDK19/CycC       | <input type="checkbox"/> EIF2AK3              | <input type="checkbox"/> GRK7             |
| <input type="checkbox"/> Bmpr1B                   | <input type="checkbox"/> CDK20/CycH       | <input type="checkbox"/> EML4 ALK             | <input type="checkbox"/> GSG2             |
| <input type="checkbox"/> BMX                      | <input type="checkbox"/> CDK20/CycT1      | <input type="checkbox"/> EML4 ALK F1174L      | <input type="checkbox"/> GSK3-alpha       |
| <input type="checkbox"/> B-RAF wt                 | <input type="checkbox"/> CHK1             | <input type="checkbox"/> EPHA1                | <input type="checkbox"/> GSK3-beta        |
| <input type="checkbox"/> B-RAF V600E              | <input type="checkbox"/> CHK2             | <input type="checkbox"/> EPHA2                | <b>H</b>                                  |
| <input type="checkbox"/> BRK                      | <input type="checkbox"/> CK1-alpha1       | <input type="checkbox"/> EPHA3                | <input type="checkbox"/> HCK              |
| <input type="checkbox"/> BRSK1                    | <input type="checkbox"/> CK1-delta        | <input type="checkbox"/> EPHA4                |                                           |
| <input type="checkbox"/> BRSK2                    | <input type="checkbox"/> CK1-epsilon      |                                               |                                           |
| <input type="checkbox"/> BTK                      | <input type="checkbox"/> CK1-gamma1       |                                               |                                           |
| <input type="checkbox"/> BUB1B                    |                                           |                                               |                                           |

# SCREENING PROJECT FORM (SPF)

## Kinases to be tested (please mark):

(list continues on following page)

- HIPK1
- HIPK2
- HIPK3
- HIPK4
- HRI

### I

- IGF1-R
- IKK-alpha
- IKK-beta
- IKK-epsilon
- INS-R
- INSR-R
- IRAK1
- IRAK4 (untagged)
- ITK

### J

- JAK1 aa583-1154wt
- JAK1 aa583-1154 S729C
- JAK2
- JAK3
- JNK1
- JNK2
- JNK3

### K

- KIT wt
- KIT V559D
- KIT V559D/V654A
- KIT V559D/T670I
- KIT V560G
- KIT V654A
- KIT T670I
- KIT D816H
- KIT D816V
- KIT A829P

### L

- LCK
- LIMK1
- LIMK2
- LRRK2 wt
- LRRK2 R1441C
- LRRK2 G2019S
- LRRK2 I2020T
- LTK
- LYN

### M

- MAP3K1
- MAP3K7/MAP3K7IP1
- MAP3K9
- MAP3K10
- MAP3K11
- MAP4K2
- MAP4K4
- MAP4K5
- MAPKAPK2
- MAPKAPK3
- MAPKAPK5
- MARK1
- MARK2
- MARK3
- MARK4
- MASTL
- MATK
- MEK1 wt
- MEK1 F53L
- MEK2
- MEK5
- MEKK2
- MEKK3
- MELK
- MERTK
- MET wt
- MET F1200I
- MET D1228H
- MET D1228N
- MET L1195V
- MET Y1230A
- MET Y1230C
- MET Y1230D
- MET Y1230H
- MET Y1235D
- MET M1250T
- MINK1
- MKK4
- MKK6 S207DT211D
- MKK7
- MKNK1
- MKNK2
- MLK4
- MST1
- MST2
- MST3
- MST4
- mTOR

- MUSK
- MYLK
- MYLK2
- MYLK3

### N

- NEK1
- NEK2
- NEK3
- NEK4
- NEK6
- NEK7
- NEK9
- NEK11
- NIK
- NLK
- NPM1 ALK
- NPM1 ALK F1174L

### P

- p38-alpha
- p38-beta
- p38-gamma
- p38-delta
- PAK1
- PAK2
- PAK3
- PAK4
- PAK6
- PAK7
- PASK
- PBK
- PDGFR-alpha wt
- PDGFR-alpha V561D
- PDGFR-alpha T674I
- PDGFR-alpha D842V
- PDGFR-beta
- PDK1
- PHKG1
- PHKG2
- PIM1
- PIM2
- PIM3
- PKA

- PKC-alpha
- PKC-beta1
- PKC-beta2
- PKC-delta
- PKC-epsilon
- PKC-eta
- PKC-gamma
- PKC-iota
- PKC-mu
- PKC-nu
- PKC-theta
- PKC-zeta
- PKC-zeta wt aa184-592 (PKM-zeta)
- PKMYT1
- PKN3
- PLK1
- PLK3
- PRK1
- PRK2
- PRKD2
- PRKG1
- PRKG2
- PRKX
- PYK2

### R

- RAF1 Y340D/Y341D (untagged)
- RET wt
- RET G691S
- RET R749T
- RET E762Q
- RET Y791F
- RET V804E
- RET V804L
- RET V804M
- RET Y806H
- RET R813Q
- RET S891A
- RET M918T
- RIPK2
- RIPK4
- RIPK5

## Kinases to be tested (please mark):

- ROCK1
- ROCK2
- RON
- ROS
- RPS6KA1
- RPS6KA2
- RPS6KA3
- RPS6KA4
- RPS6KA5
- RPS6KA6

### S

- S6K
- S6K-beta
- SAK
- SGK1
- SGK2
- SGK3
- SIK1
- SIK2
- SIK3
- SLK
- SNARK
- SNK
- SRC (GST-HIS-tag)
- SRMS
- SRPK1
- SRPK2
- STK17A
- STK23
- STK25
- STK33
- STK39
- SYK aa1-635

### T

- TAOK2
- TAOK3
- TBK1
- TEC
- TGFB-R1
- TGFB-R2
- TIE2 wt
- TIE2 R849W
- TIE2 Y897S
- TIE2 Y1108F

- TLK1
- TLK2
- TNK1
- TRK-A wt
- TRK-A G667C
- TRK-B
- TRK-C
- TSF1
- TSK2
- TSSK1
- TTBK1
- TTBK2
- TTK
- TXK
- TYK2
- TYRO3

### U

- ULK2

### V

- VEGF-R1
- VEGF-R2
- VEGF-R3
- VRK1
- VRK2

### W

- WEE1
- WNK1
- WNK2
- WNK3

### Y

- YES

### Z

- ZAK
- ZAP70

## II. Lipid Kinases

### PI4K Class I

- PIK3CA/PIK3R1
- PIK3CB wt/PIK3R1
- PIK3CB E633K/PIK3R1
- PIK3CB E1051K/PIK3R1
- PIK3CB D1067A/PIK3R1
- PIK3CB D1067V/PIK3R1
- PIK3CB D1067Y/PIK3R1
- PIK3CB L1049R/PIK3R1
- PIK3CD/PIK3R1
- PIK3CG

### PI4K Class II

- PIK3C2A
- PIK3C2B
- PIK3C2G
- PI4K2A
- PI4K2B

### PI4K Class III

- PIK3C3

### PI4K Class IV

- PI4KB

### PIP5K1

- PIP5K1A
- PIP5K1B
- PIP5K1C

# SCREENING PROJECT FORM (SPF)

## Check Lists

### A) General check list:

Compound Submission Form (vials) or file with compound coordinates (plates) mailed to f.totzke@proqinase.com?

A copy of this Screening Project Form (dated & signed) included into package?

### B) Check list for dry ice plate shipment:

Plate(s) properly sealed and lidded (see guidelines)?

If more than one plate: Plates sealed & stacked, upper-most plate(s) lidded, stack(s) wrapped with rubber band and stack(s) enclosed in plastic bag(s)?

Sufficient dry ice included (at least 5 kg/prospective day of transport)?

## Signature

Date

Signature

## Shipping Address

**Dr. Frank Totzke**

**Biochemical Screening Systems**

**ProQinase GmbH**

**Breisacher Strasse 117**

**79106 Freiburg**

**Germany**

**phone: +49-761-769996-1723**

**fax: +49-761-769996-1781**

**mail: f.totzke@proqinase.com**