

Please fill in the parts shown in green background



NanoBRET™ Target Engagement Intracellular Kinase Cell-Based Assay Services

Please read the "Important Notice" stated below thoroughly before proceeding.

Rev. 2018_11

Date : Enter mm/dd/yy

Customer Information [Highlighted fields are required.]

Customer Name
Institution/Company
Department
Street Address 1
Street Address 2
City
State/Zip
Country
Tel Number
Fax Number
Email Address

Dr. John Carna

Japan

Enter Your Information

Attention

Cells highlighted in blue require information and will be colored according to the information you type.
Please do not delete any cells in this spreadsheet.

Assay Information

IC₅₀ Determination

Eight (8) concentrations tested in half-log increments

Technical Replicates

☐ n=1
☐ n=2
☐ Others

Choose the number of your requested replication

Compound Information [Highlighted fields are required.]

Compound State

☒ Solution

☐ Solid

No.	Compound Name	Conc. (mM)	Volume (μL)	Purity (%)	Storage Temperature	Please fill in the Highest Test Concentration				
1	Enter information on one (1) compound per line				Freeze	Conc 1. (μM)				
2						10	3, 1	0.3, 01	0.03, 001	0.003
3							0, 0	0, 0	0, 0	0
4							0, 0	0, 0	0, 0	0
5							0, 0	0, 0	0, 0	0
6							0, 0	0, 0	0, 0	0
7							0, 0	0, 0	0, 0	0
8							0, 0	0, 0	0, 0	0
9							0, 0	0, 0	0, 0	0
10							0, 0	0, 0	0, 0	0
11							0, 0	0, 0	0, 0	0
12							0, 0	0, 0	0, 0	0
13							0, 0	0, 0	0, 0	0
14							0, 0	0, 0	0, 0	0
15							0, 0	0, 0	0, 0	0

Note: Your compound provided for this service will be disposed 12 weeks after completion of study. For other options, please consult us.

Target Kinase Selection

NanoBRET™ TE Intracellular Kinase Assay Panel

(Bold letters: Off-the-Shelf targets, Small letters: Custom targets)

Total 0 target(s) selected

(Off-the-Shelf 0 targets/Custom 0 targets)

- ☐ AAK1

☐ ABL1

☐ ABL1 [E255K]

☐ ABL1 [F317I]

☐ ABL1 [F317L]

☐ ABL1 [H396P]

☐ ABL1 [M351T]

☐ ABL1 [Q252H]

☐ ABL1 [Y253F]

☐ ABL2

☐ ACVR1

☐ ACVR1B

☐ ACVR1L1

☐ ADK

☐ AKT2

☐ ALK [L1196M]

☐ AMHR2

☐ AURKA

☐ AURKB

☐ AURKC

☐ AXL

☐ BLK

☐ BMP2K

☐ BMPR1A

☐ BMX

☐ BRSK1

☐ BRSK2

☐ BTK

☐ CAMK2A

☐ CAMK2D

☐ CAMK2G

☐ CDK1/Cyc B1

☐ CDK1/Cyc B2

☐ CDK1/Cyc E1

☐ CDK1/Cyc K

☐ CDK10/Cyc L2

☐ CDK11A/Cyc K

☐ CDK11A/Cyc L2

☐ CDK14/Cyc Y

☐ CDK15/Cyc Y

☐ CDK16/Cyc Y

☐ NTRC Oncolines™

☐ CSNK1D

☐ CSNK1E

☐ CSNK1

☐ CSNK2A1

☐ CSNK2A2

☐ DAPK2

☐ DCLK1

☐ DCLK3

☐ DDR1

☐ DDR2

☐ DYRK1A

☐ DYRK1B

☐ EIF2AK4k2

☐ EPHA1

☐ EPHA2

☐ EPHA4

☐ EPHA5

☐ EPHA6

☐ EPHA7

☐ EPHA8

☐ EPHB1

☐ EPHB2

☐ EPHB3

☐ EPHB4

☐ ERN1

☐ ERN2

☐ FER

☐ FES

☐ FGFR1

☐ FGFR2

☐ FGFR3

☐ FGFR4

☐ FGR

☐ FLT1

☐ FLT3

☐ FLT3 [D835H]

☐ FLT3 [D835V]

☐ FLT3 [D835Y]

☐ FLT3 [K663Q]

☐ FLT3 [N841I]

☐ FLT3 [R834Q]

☐ LRRK2

☐ LRRK2 [G2019S]

☐ LTK

☐ LYN

☐ MAP3K10

☐ MAP3K11

☐ MAP3K12

☐ MAP3K19

☐ MAP3K2

☐ MAP3K21

☐ MAP3K3

☐ MAP3K4

☐ MAP3K9

☐ MAP4K1

☐ MAP4K2

☐ MAP4K3

☐ MAP4K5

☐ MAPK1

☐ MAPK11

☐ MAPK13

☐ MAPK14

☐ MAPK3

☐ MAPK4

☐ MAPK6

☐ MAPK8

☐ MAPK9

☐ MARK2

☐ MARK3

☐ MARK4

☐ MAST4

☐ MELK

☐ MET

☐ MET [M1250T]

☐ MET [Y1235D]

☐ MKNK1

☐ MKNK2

☐ MLTK

☐ MOK

☐ MUSK

☐ MYLK2

☐ MYLK3

☐ PIKFYVE

☐ PIP4K2C

☐ PIP5K1B

☐ PKMYT1

☐ PLK2

☐ PLK3

☐ PLK4

☐ PRKAA2

☐ PRKACA

☐ PRKACB

☐ PRKCE

☐ PRKG2

☐ PRKX

☐ PTK2

☐ PTK2B

☐ PTK6

☐ RET

☐ RET [M918T]

☐ RET [V804L]

☐ RET [V804M]

☐ R1OK2

☐ RIPK1

☐ RIPK2

☐ RON

☐ RPS6KA1

☐ RPS6KA2

☐ RPS6KA3

☐ RPS6KA4

☐ RPS6KA6

☐ SBK3

☐ SGK1

☐ SIK1

☐ SIK3

☐ SLK

☐ SNF1LK2

☐ SNRK

☐ SRC

☐ SRMS

☐ STK10

☐ STK16

☐ STK3

Instruction for Shipping Compound <Important>

Minimum required volume for your study: 50 uL

Required compound volume is shown here

Number of targets	Required compound volume
1-8	50 uL
8-32	100 uL
33-80	200 uL
81-160	300 uL
161-295	500 uL

Please ship your compound(s) to

c/o Carna Biosciences, Inc.
BMA3F 1-5-5 Minatojima-Minamimachi
Chuo, Kobe 650-0047 Japan

E-mail: info@carnabio.com

Please double check before you ship...

- * **Minimum required volume**
- * **1,000 X highest test conc. dissolved in 100% DMSO**
- * **Labeled and tightly sealed**
- * **Shipment arranged with enough dry ice and assigned date**

Additional Information**If any, please fill in****Important Notice****Turnaround**

Turnaround time is 3 weeks after determining test concentration(s) of your compound(s).

Intellectual Property and Deliverables

All work products resulting from this services shall be the sole and exclusive property of the client.

Ship-To Address

Ship test sample with the concentration and volume auto-calculated in the section of "Compound Information" above via appropriate method.

BMA 3F, 1-5-5, Minatojima-Minamimachi, Chuo-ku,
Kobe JAPAN 650-0047
TEL: 078-302-7091 / FAX: 078-302-7086

To prevent leakage and/or damage, please package your test sample securely.
Carna takes no responsibility for any damage or loss caused in transportation.
You will be notified via E-mail or FAX when your sample arrived at our facility.

Residuals

Residuals will be disposed after 12 weeks from completion of the study.
Should you have any questions or concerns, please feel free to contact us.