



ResidenceTimer™ Target Residence Time Measurements

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

Rev. 2014_12

本サービスの申込用紙は「英語版」のみです。予めご了承下さい。

Date : 申込日

Customer Information [Highlighted fields are required.]

Customer Name	カルナ太郎
Institution/Company	
Department	
Street Address 1	
Street Address 2	
City	
State/Zip	
Country	Japan
Tel Number	
Fax Number	
Email Address	

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.

お客様情報をご記入ください。
色付きのセルには必ずご記入ください。

Assay Information

Affinity Determination - Binding and Dissociation Kinetics
Five (5) concentrations tested in square root (10) dilutions

Replicates

- ☐ Single Measurement
☐ Duplicate ご希望のn数をチェックください。
☐ Triplicate

Compound Information [Highlighted fields are required.]

No.	Name of Compound	Storage Temp	Highest Test Conc.			Required Stock Conc.	Required Volume
			10 uM	1 uM	100 nM		
1	CBS-001		●	●	●	10,000 uM	100 uL
2			○	○	○	10,000 uM	100 uL
3			○	○	○	10,000 uM	100 uL
4			○	○	○	10,000 uM	100 uL
5			○	○	○		
6			○	○	○		
7			○	○	○		
8			○	○	○		
9			○	○	○		
10			○	○	○	10,000 uM	100 uL
11			○	○	○	10,000 uM	100 uL
12			○	○	○	10,000 uM	100 uL
13			○	○	○	10,000 uM	100 uL
14			○	○	○	10,000 uM	100 uL
15			○	○	○	10,000 uM	100 uL

ご希望のトップ試験濃度を選択してください。
お送りいただく被験物質の濃度、ならびに最低必要量が自動で表示されます。

Note: Your compound provided for this service is NOT returnable and will be disposed three (3) month after completion of study.

Target Kinases

試験したいキナーゼにチェックマークを付けてください。選択されるとキナーゼ名が赤色に変わります。

NTRC's Validated Kinases for ResidenceTimer™

checked or RED colored kinases are selected.

☐ ResidenceTimer™
Full Panel
(42 kinases)

フルパネルをご希望の場合は上記ボックスに
チェックマークを付けてください。

Biotinylated Kinases

☒ ABL (ABL1)

- ☐ AKT1
☐ AKT2
☐ ALK
☐ ALK[L1196M]
☐ AurA (AURKA)
☐ AXL
☐ BMX
☐ BTK
☐ CHK1 (CHEK1)
☐ DDR2
☐ EGFR
☐ FAK (PTK2)
☐ FGFR1
☐ FGFR3
☐ FGFR4
☐ FLT3
☐ FMS (CSF1R)
☐ FYN [isoform a]
☐ IGF1R
☐ IRAK4
☐ ITK
☐ KDR (VEGFR2)
☐ LCK
☐ LYNa
☐ MAP2K1

Kinase Domain

- Full length
Catalytic domain
Catalytic domain
Cytoplasmic domain
Cytoplasmic domain
Full length
Cytoplasmic domain
Full length
Full length
Full length
Full length
Cytoplasmic domain
Cytoplasmic domain
Truncated human
Cytoplasmic domain
Cytoplasmic domain
Cytoplasmic domain
Cytoplasmic domain
Full length
Cytoplasmic domain
Full length
Full length
Cytoplasmic domain
Full length
Full length
Full length

Biotinylated Kinases

- ☐ MAP3K5
☐ MET
☐ p38α (MAPK14)
☐ p38α (MAPK14)
☐ PAK4
☐ PIK3CA/PIK3R1
☐ PKN1
☐ PYK2 (PTK2B)
☐ RET
☐ ROCK1
☐ SRC
☐ TIE2
☐ TNIK
☐ TRKA (NTRK1)
☐ TRKB (NTRK2)
☐ TRKC (NTRK3)

Kinase Domain

- Catalytic domain
Cytoplasmic domain
Truncated human
Truncated human
Full length
Full length
Full length
Full length
Cytoplasmic domain
Catalytic domain
Full length
Cytoplasmic domain
Catalytic domain
Cytoplasmic domain
Cytoplasmic domain
Catalytic domain

Total 1 Kinases Selected

Instruction for Shipping Compound <Important>

Minimum required volume for your study 100 uL

Number of kinases selected	Minimum required volume
1-42	100 uL

Please double check before you ship...

- * Minimum required volume
- * Labeled and tightly sealed
- * Shipment arranged with enough dry ice and assigned date

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

Additional Information

Important Notice (ご注意事項)

納期

Turnaround

納期は試験開始後、約3週間です。(SynergyFinder™は含まず。)
Turnaround time is approximately 3 weeks, excluding SynergyFinder™.

成果の帰属

Intellectual Property and Deliverables

本サービスにおける成果物(データ)は試験委託者に帰属し、試験委託者はこれを自由に利用または公表することができます。
また、成果物および成果物に基づいて得られる知的財産権は試験委託者に帰属します。
All work products resulting from this services shall be the sole and exclusive property of the client.

試験物質の送付先

Ship-To Address

被験物質は「Compound Information」欄に自動表示された指定の濃度、容量を適切な輸送方法にて下記宛にお送り下さい。
Ship test sample with the concentration and volume auto-calculated in the section of "Compound Information" above via appropriate method.

〒650-0047 神戸市中央区港島南町1丁目5番5号 BMA 3F
カルナバイオサイエンス株式会社 営業部 宛
TEL: 078-302-7091 / FAX: 078-302-7086

輸送中にこぼれたり、容器が破損したりしないよう梱包には十分にご注意ください。
輸送中における被験物質のトラブルに関して弊社では責任を負いかねますのでご了承ください。
被験物質を受領後、弊社よりメール、またはファクシミリにて受領確認のご連絡を差し上げます。
お預かりした被験物質は厳重に梱包の上、NTRC社に発送いたします。
To prevent leak and/or damage, please pack your test sample securely.
Carna takes no responsibility for any damage or loss caused in transit.
You will be notified via E-mail or FAX when your sample arrived at our facility.
Your test sample will be forwarded to NTRC with care.

残余被験物質

Residuals

お預かりしました被験物質は、原則、最終報告書を提出いたしました日の3ヵ月後に廃棄させていただきます。
その他、ご不明な点・ご要望等ございましたら、ご遠慮なくお申し付けください。
Residuals will be disposed after three (3) months from completion of the study.
Should you have any questions or concerns, please feel free to contact us.