

Biotinylated Kinases

65
種類

裏面リストを
ご覧ください

表面プラズモン共鳴(SPR)やバイオレイヤー干渉法(BLI)といった物質間の相互作用を評価する系(解析機器)でご利用いただけるビオチン化キナーゼ蛋白質を販売しております。これらの解析機器で低分子化合物の創薬研究・開発を行なう際、リガンドとする標的分子蛋白質を活性、構造を保持したままセンサー表面に固定化(固相化)することはとても難しいことです。当社はこれまで300を超えるキナーゼ種を自社製造してきた上で培ったノウハウを元に下記の特長をもつビオチン化キナーゼ蛋白質の開発成功に至り、販売する運びとなりました。

ビオチン化キナーゼ蛋白質

このような特長や利点があります

- 一つのキナーゼ分子に一つのビオチン分子をラベリング
- 特異的部位をビオチン化
- バキュロウイルス発現システムにより蛋白質を取得
- 購入後のラベリングが不要で取扱いが簡単 → 時間、コスト削減に貢献



ビオチン化キナーゼ蛋白質(FLT3)の使用例

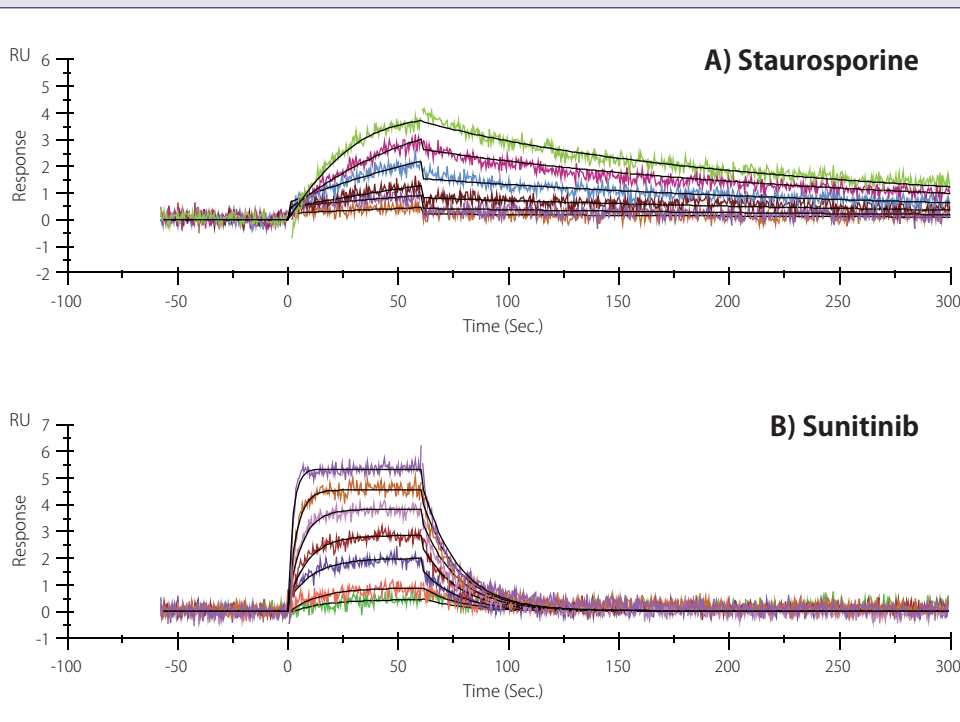


図 1. Representative SPR sensorgrams of the binding between biotinylated-FLT3 and kinase inhibitors. A) Staurosporine, and B) Sunitinib.

製造ステップ

クローニング、培養



精製

- 未処理

将来的に以下の処理オプション(有料)をお客様に選択いただけるよう現在検討を進めています

- ATP 処理
- 脱リン酸化処理

※ 詳細はお問合せ下さい



品質チェック



入庫

販売単位: 10µg/100µg

SPR(表面プラズモン共鳴)やBLI(バイオレイヤー干渉法)だけでなく
様々な用途、解析機器でご利用いただける製品です

カルナバイオサイエンス株式会社

〒650-0047 神戸市中央区港島南町1-5-5 BMA3F

TEL: 078-302-7091 (営業部直通) / FAX: 078-302-7086

E-mail: info@carnabio.com

Biotinylated Kinases from Carina Biosciences, Inc.

Cat No.	Name of Product	Length	Amino Acid	Accession Number	Tag	M.W.	Expression System
08-401-20N	BTN - ABL(ABL1)	Full-length	2-1130(end)	NP_005148.2	N-terminal DYKDDDDK	126 kDa	Insect (<i>sf21</i>)
09-432-20N	BTN - ACTR2B(ACVR2B)	Cytoplasmic domain	161-512(end)	NP_001097.2	N-terminal DYKDDDDK	43 kDa	Insect (<i>sf21</i>)
01-401-20N	BTN - AKT1	Catalytic domain	104-480(end)	NP_005154.1	N-terminal DYKDDDDK	47 kDa	Insect (<i>sf21</i>)
01-402-20N	BTN - AKT2	Catalytic domain	120-481(end)	NP_001617.1	N-terminal DYKDDDDK	45 kDa	Insect (<i>sf21</i>)
08-405-20N	BTN - ALK	Cytoplasmic domain	1058-1620(end)	BAG10812.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>sf21</i>)
08-429-20N	BTN - ALK[L1196M]	Cytoplasmic domain	1058-1620(end)	BAG10812.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>sf21</i>)
05-401-20N	BTN - AurA(AURKA)	Full-length	1-403(end)	NP_940835.1	N-terminal DYKDDDDK	50 kDa	Insect (<i>sf21</i>)
05-402-21N	BTN - AurB(AURKB)/INCENP	Full-length	1-344	Q96GD4	N-terminal DYKDDDDK	43 kDa	Insect (<i>sf21</i>)
08-407-20N	BTN - AXL	Cytoplasmic domain	464-885(end)	NP_001690.2	N-terminal DYKDDDDK	51 kDa	Insect (<i>sf21</i>)
08-479-20N	BTN - BMX	Full-length	1-675(end)	NP_001712.1	N-terminal DYKDDDDK	82 kDa	Insect (<i>sf21</i>)
08-480-20N	BTN - BTK	Full-length	2-659(end)	NP_000052.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>sf21</i>)
02-417-20N	BTN - CHK1(CHEK1)	Full-length	1-476(end)	NP_001265.1	N-terminal DYKDDDDK	58 kDa	Insect (<i>sf21</i>)
08-414-20N	BTN - DDR2	Cytoplasmic domain	422-855(end)	NP_006173.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)
09-411-20N	BTN - DLK(MAP3K12)	Catalytic domain	1-520	NP_006292.3	N-terminal DYKDDDDK	63 kDa	Insect (<i>sf21</i>)
08-415-20N	BTN - EGFR	Cytoplasmic domain	669-1210(end)	NP_005219.2	N-terminal DYKDDDDK	66 kDa	Insect (<i>sf21</i>)
08-410-20N	BTN - EGFR[T790M/L858R]	Cytoplasmic domain	669-1210(end)	NP_005219.2	N-terminal DYKDDDDK	66 kDa	Insect (<i>sf21</i>)
08-437-20N	BTN - FAK(PTK2)	Truncated protein	376-1052(end)	NP_722560.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>sf21</i>)
08-435-20N	BTN - FGFR1	Cytoplasmic domain	396-820(end)	NP_056934.2	N-terminal DYKDDDDK	52 kDa	Insect (<i>sf21</i>)
08-433-20N	BTN - FGFR3	Cytoplasmic domain	436-806(end)	NP_000133.1	N-terminal DYKDDDDK	52 kDa	Insect (<i>sf21</i>)
08-436-20N	BTN - FGFR4	Cytoplasmic domain	460-802(end)	NP_002002.3	N-terminal DYKDDDDK	42 kDa	Insect (<i>sf21</i>)
08-454-20N	BTN - FLT3	Cytoplasmic domain	564-993(end)	NP_004110.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)
08-455-20N	BTN - FMS(CSF1R)	Cytoplasmic domain	538-972(end)	NP_005202.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)
08-468-20N	BTN - FYN[isoform a]	Full-length	1-537(end)	NP_002028.1	N-terminal DYKDDDDK	65 kDa	Insect (<i>sf21</i>)
08-441-20N	BTN - IGF1R	Cytoplasmic domain	959-1367(end)	NP_000866.1	N-terminal DYKDDDDK	50 kDa	Insect (<i>sf21</i>)
05-484-20N	BTN - IKKβ(IKKB)	Truncated protein	1-662	NP_001547.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>sf21</i>)
05-412-20N	BTN - IKKα(CHUK)	Full-length	1-745(end)	NP_001269.3	N-terminal DYKDDDDK	88 kDa	Insect (<i>sf21</i>)
05-412-11-20N	BTN - IKKα(CHUK)[inactive mutant]	Full-length	1-745(end)	NP_001269.3	N-terminal DYKDDDDK	88 kDa	Insect (<i>sf21</i>)
09-445-20N	BTN - IRAK4	Full-length	1-460(end)	NP_057207.2	N-terminal DYKDDDDK	55 kDa	Insect (<i>sf21</i>)
08-481-20N	BTN - ITK	Full-length	2-620(end)	NP_005537.3	N-terminal DYKDDDDK	75 kDa	Insect (<i>sf21</i>)
08-444-20N	BTN - JAK1	Catalytic domain	850-1154(end)	NP_002218.2	N-terminal DYKDDDDK	39 kDa	Insect (<i>sf21</i>)
08-445-20N	BTN - JAK2	Catalytic domain	826-1132(end)	NP_004963.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>sf21</i>)
08-491-20N	BTN - KDR(VEGFR2)	Cytoplasmic domain	790-1356(end)	NP_002244.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>sf21</i>)
08-456-21N	BTN - KIT	Cytoplasmic domain	544-976(end)	NP_000213	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)
08-470-20N	BTN - LCK	Full-length	1-509(end)	NP_005347.2	N-terminal DYKDDDDK	62 kDa	Insect (<i>sf21</i>)
08-471-20N	BTN - LYNa	Full-length	1-512(end)	NP_002341.1	N-terminal DYKDDDDK	63 kDa	Insect (<i>sf21</i>)
07-441-10-20N	BTN - MAP2K1[inactive]	Full-length	1-393(end)	NP_002746.1	N-terminal DYKDDDDK	47 kDa	Insect (<i>sf21</i>)
07-407-20N	BTN - MAP3K5	Catalytic domain	654-971	NP_005914.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>sf21</i>)
08-451-20N	BTN - MET	Cytoplasmic domain	956-1390(end)	NP_000236.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)
02-445-20N	BTN - MNK1(MKNK1)	Full-length	1-424(end)	BAA19885.1	N-terminal DYKDDDDK	51 kDa	Insect (<i>sf21</i>)
05-433-20N	BTN - NEK9	Truncated protein	1-346, 733-979(end)	NP_149107.4	N-terminal DYKDDDDK	70 kDa	Insect (<i>sf21</i>)
04-452-20N	BTN - p38α(MAPK14)	Truncated protein	9-352, 1-334(end)	NP_620581.1	N-terminal DYKDDDDK	43 kDa	Insect (<i>sf21</i>)
04-452-10-20N	BTN - p38α(MAPK14)[inactive]	Truncated protein	9-352	NP_620581.1	N-terminal DYKDDDDK	43 kDa	Insect (<i>sf21</i>)
07-426-20N	BTN - PAK4	Full-length	1-591(end)	NP_005875.1	N-terminal DYKDDDDK	68 kDa	Insect (<i>sf21</i>)
11-401-20N	BTN - PIK3CA/PIK3R1	Full-length	1-1068(end)	NP_006209.2	N-terminal DYKDDDDK	128 kDa	Insect (<i>sf21</i>)
01-438-20N	BTN - PKCη(PRKCH)	Full-length	1-683(end)	NP_006246	N-terminal DYKDDDDK	82 kDa	Insect (<i>sf21</i>)
01-444-20N	BTN - PKN1	Full-length	1-942(end)	NP_002732.3	N-terminal DYKDDDDK	108 kDa	Insect (<i>sf21</i>)
05-457-20N	BTN - PLK1	Full-length	1-603(end)	NP_005021.2	N-terminal DYKDDDDK	72 kDa	Insect (<i>sf21</i>)
08-438-20N	BTN - PYK2(PTK2B)	Full-length	1-967(end)	NP_775267.1	N-terminal DYKDDDDK	115 kDa	Insect (<i>sf21</i>)
08-459-20N	BTN - RET	Cytoplasmic domain	658-1114(end)	NP_066124.1	N-terminal DYKDDDDK	56 kDa	Insect (<i>sf21</i>)
01-409-20N	BTN - ROCK1	Catalytic domain	1-477	NP_005397.1	N-terminal DYKDDDDK	59 kDa	Insect (<i>sf21</i>)
11-405-20N	BTN - SPHK1	Full-length	1-384(end)	NP_001136074.1	N-terminal DYKDDDDK	46 kDa	Insect (<i>sf21</i>)
08-473-20N	BTN - SRC	Full-length	1-536(end)	NP_005408.1	N-terminal DYKDDDDK	64 kDa	Insect (<i>sf21</i>)
04-460-20N	BTN - SRPK1	Full-length	1-654(end)	CAI20544.1	N-terminal DYKDDDDK	78 kDa	Insect (<i>sf21</i>)
11-406-20N	BTN - SRPK2	Full-length	1-618(end)	NP_001191089.1	N-terminal DYKDDDDK	69 kDa	Insect (<i>sf21</i>)
08-476-20N	BTN - SYK	Full-length	1-635(end)	NP_003168.2	N-terminal DYKDDDDK	76 kDa	Insect (<i>sf21</i>)
09-419-21C	BTN - TAK1-TAB1(MAP3K7)	Fused protein	1-303	NP_663304.1	C-terminal DYKDDDDK	45 kDa	Insect (<i>sf21</i>)
09-441-20N	BTN - TGFβR1(TGFB1)	Catalytic domain	200-503(end)	NP_004603.1	N-terminal DYKDDDDK	38 kDa	Insect (<i>sf21</i>)
08-485-20N	BTN - TIE2	Cytoplasmic domain	771-1124(end)	NP_000450.1	N-terminal DYKDDDDK	44 kDa	Insect (<i>sf21</i>)
07-438-20N	BTN - TNIK	Catalytic domain	1-314	NP_055843.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>sf21</i>)
08-486-20N	BTN - TRKA(NTRK1)	Cytoplasmic domain	436-790(end)	NP_001012331.1	N-terminal DYKDDDDK	44 kDa	Insect (<i>sf21</i>)
08-487-20N	BTN - TRKB(NTRK2)	Cytoplasmic domain	456-822(end)	NP_001018074.1	N-terminal DYKDDDDK	45 kDa	Insect (<i>sf21</i>)
08-497-20N	BTN - TRKC(NTRK3)	Catalytic domain	456-825(end)	NP_002521.2	N-terminal DYKDDDDK	46 kDa	Insect (<i>sf21</i>)
08-447-20N	BTN - TYK2	Catalytic domain	871-1187(end)	NP_003322.3	N-terminal DYKDDDDK	40 kDa	Insect (<i>sf21</i>)
05-479-20N	BTN - WNK1	Catalytic domain	1-491	NP_061852.1	N-terminal DYKDDDDK	59 kDa	Insect (<i>sf21</i>)
05-482-20N	BTN - WNK4	Catalytic domain	1-444	NP_115763.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>sf21</i>)

Updated: 2015/4/14

お知らせ：掲載の内容は予告なく変更する場合があります。予めご了承下さい。