

Biotinylated Kinases

Biotinylated Kinases

Biotinylated kinases for the study of kinase-molecule interactions are now available from Carna Biosciences.

The immobilization of target proteins onto sensor surfaces, while maintaining both structure and activity, has been problematic and rate-limiting to the acquisition of meaningful SPR and/or BLI data in small molecule drug discovery. Carna Biosciences has developed proprietary technology to produce biotinylated kinase products to help overcome these challenges. A few of this new and expanding product line are featured below.

Biotinylated Kinases... The Carna Advantage

- ➡ **Biotinylation at a single site outside the catalytic domains;**
- ➡ **Enzymatic activity and structure are retained;**
- ➡ **Baculovirus expression system (insect cells) used for production;**
- ➡ **Carna's biotinylated kinases are highly purified by affinity chromatographic procedures.**

Released October 2010

The table below lists biotinylated kinases just released! Additional biotinylated kinases under development are listed on the reverse side. Please contact us for the availability of products not yet listed.

A sample product sheet is available at our website, www.carnabio.com or contact your area representative.

Product Name		Catalog No.	List Price - 10µg	List Price - 100µg
BTN - ABL(ABL1)	Cytoplasmic Tyrosine	08-401-20N	US\$ 1,400	US\$ 8,000
BTN - AKT1	Serine/Threonine	01-401-20N	US\$ 1,400	US\$ 8,000
BTN - ALK	Receptor Tyrosine	08-405-20N	US\$ 1,400	US\$ 8,000
BTN - AXL	Receptor Tyrosine	08-407-20N	US\$ 1,400	US\$ 8,000
BTN - BTK	Cytoplasmic Tyrosine	08-480-20N	US\$ 1,400	US\$ 8,000
BTN - DLK(MAP3K12)	Serine/Threonine	09-411-20N	US\$ 1,400	US\$ 8,000
BTN - EGFR	Receptor Tyrosine	08-415-20N	US\$ 1,400	US\$ 8,000
BTN - FGFR1	Receptor Tyrosine	08-435-20N	US\$ 1,400	US\$ 8,000
BTN - FMS(CSF1R)	Receptor Tyrosine	08-455-20N	US\$ 1,400	US\$ 8,000
BTN - HER2(ERBB2)	Receptor Tyrosine	08-416-20N	US\$ 1,400	US\$ 8,000
BTN - IGF1R	Receptor Tyrosine	08-441-20N	US\$ 1,400	US\$ 8,000
BTN - IKK α (CHUK)	Serine/Threonine	05-412-20N	US\$ 1,400	US\$ 8,000
BTN - JAK1	Cytoplasmic Tyrosine	08-444-20N	US\$ 1,400	US\$ 8,000
BTN - JAK2	Cytoplasmic Tyrosine	08-445-20N	US\$ 1,400	US\$ 8,000
BTN - KDR	Receptor Tyrosine	08-491-20N	US\$ 1,400	US\$ 8,000
BTN - MAP2K1	Serine/Threonine	07-441-20N	US\$ 1,400	US\$ 8,000
BTN - MET	Receptor Tyrosine	08-451-20N	US\$ 1,400	US\$ 8,000
BTN - PIK3CA/PIK3R1	Lipid	11-401-20N	US\$ 1,400	US\$ 8,000
BTN - PLK1	Serine/Threonine	05-457-20N	US\$ 1,400	US\$ 8,000
BTN - PLK4	Serine/Threonine	05-460-20N	US\$ 1,400	US\$ 8,000
BTN - SYK	Cytoplasmic Tyrosine	08-476-20N	US\$ 1,400	US\$ 8,000
BTN - TGF β R1(TGFBR1)	Serine/Threonine	09-441-20N	US\$ 1,400	US\$ 8,000
BTN - TNIK	Serine/Threonine	07-438-20N	US\$ 1,400	US\$ 8,000

**Introductory
Promotion**

Special pricing is
now available for all
currently available
biotinylated kinases!

(Please see reverse side)

2010.11.24

NEW

2010.11.24

NEW

Note: The contents of this list are subject to change without notice.

Biotinylated Kinases

Production Steps

Cloning, Culturing Cells

Purification

■ **Standard Production**

■ **ATP-treated and**

■ **Phosphatase Treated Kinases**

will be offered on a fee-for-service basis in the future. Please contact us for more detail.

Quality Control

Off The Shelf

Vial Sizes : 10 μ g / 100 μ g

Example Data

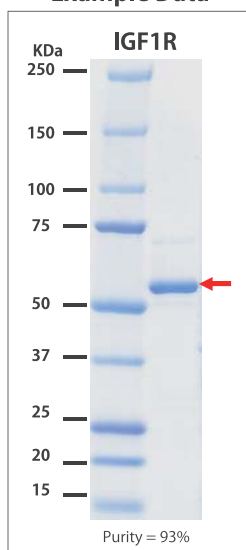


Fig1. IGF1R analysis by SDS-PAGE

Biotinylated Kinase - Example of Use

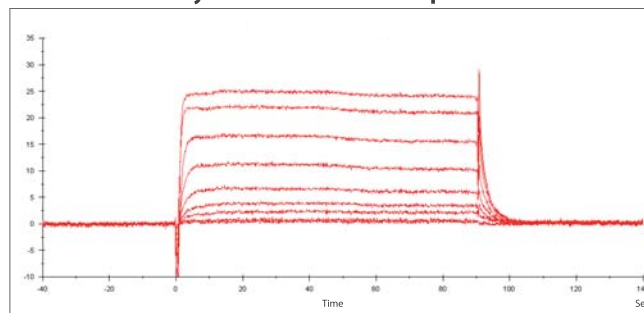


Fig2. IGF1R biotinylated kinase vs. Staurosporine binding detected using Surface Plasmon Resonance (SPR)

Under Development

Please contact us for availability of kinases not listed below.

Product Name		Catalog No.	List Price - 10 μ g	List Price - 100 μ g
BTN - BRAF	Serine/Threonine	TBD	TBD	TBD
BTN - COT(MAP3K8)	Serine/Threonine	TBD	TBD	TBD
BTN - FLT3	Receptor Tyrosine	TBD	TBD	TBD
BTN - JAK3	Cytoplasmic Tyrosine	TBD	TBD	TBD
BTN - MAP3K14	Serine/Threonine	TBD	TBD	TBD
BTN - RAF1	Serine/Threonine	TBD	TBD	TBD
BTN - TRKA	Receptor Tyrosine	TBD	TBD	TBD
BTN - TRKB	Receptor Tyrosine	TBD	TBD	TBD
BTN - TRKC	Receptor Tyrosine	TBD	TBD	TBD
BTN - TYK2	Cytoplasmic Tyrosine	TBD	TBD	TBD

Note: The contents of this list are subject to change without notice.

Introductory Promotion

Special pricing is now available for all currently available biotinylated kinases!

Promotional Pricing on **Biotinylated Kinases**

Limited Time Offer! All orders must be received by December 17, 2010 to receive our special pricing.

		List Price	Promotional Price
Discovery Pack	10 μ g	USD1,400	USD 950
Development Pack	100 μ g	USD 8,000	USD 5,500



Contact us for availability or further information
Call today 1 888 645 1233!

Carna's wholly-owned subsidiary

CarnaBio USA, Inc.

209 West Central Street, Suite 127, Natick, MA 01760 USA

TEL: +1-508-650-1244 / Toll-Free: +1-888-645-1233

FAX: +1-508-650-1722 / E-Mail: orders@carnabio.com

www.carnabio.com