

Crystallography

New Tools and Crystallography Services for Structure Based Drug Discovery

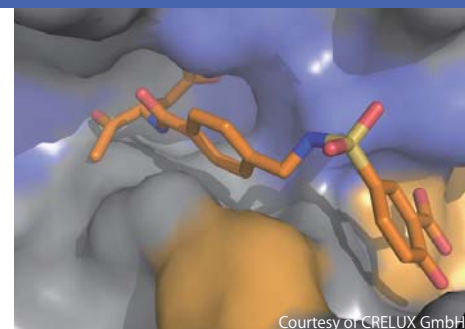


Carna now offers Co-Crystallography Services & Crystal Grade Proteins through collaboration with CRELUX GmbH. We have expanded our extensive offering of products and services to include support for Structure-based Drug Discovery. Opt for Co-crystallography Services to visualize your drug candidates with the targeted enzyme, or grow-your-own crystals using our Crystal Grade Proteins.

Targets listed on reverse side

CRELUX's Co-Crystallography Services "XPRESS"

As the process of Structure Based Drug Discovery is very important, we offer the CRELUX's crystallography service called **XPRESS** for a variety of kinases as listed on reverse side. Our collaboration project plans display a modular set up including free to choose exit points. You will profit from low fixed costs and predominantly success-based payment schedules. We are also going to offer a custom-crystallize service including DNA cloning, expression, and co-crystallization.



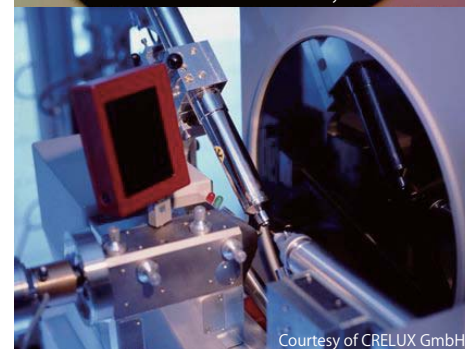
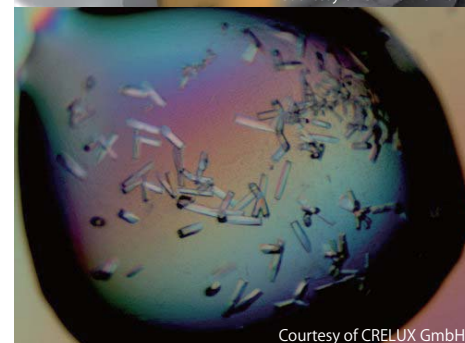
Crystal Grade Proteins (Kinase and Non-Kinase Targets)

In addition to the Co-Crystallography Service, most of Off The Shelf proteins^(*1) shown on the list can be offered as Crystal Grade Proteins^(*2) for a fast, reliable, and affordable procedure of co-crystallization performed by customer-self. Expression, purification, and crystallization for **XPRESS** targets have been established by CRELUX.

Also, crystal grade proteins developed by Carna Biosciences^(*3) solely are available. Crystallization is confirmed on all crystal grade proteins from CRELUX and Carna before dispatch. A protocol^(*4) including the condition comes with products.



Sample - CRELUX's Data Sheet



(*1) Some proteins may require compounds at the time of expression, therefore, not available for individual purchase.

(*2) Proteins may not be active, crystallization is prioritized.

(*3) Crystal grade proteins produced by Carna Biosciences has activity.

(*4) This is to check the quality of product.

For more details, please visit our website.

Carna Biosciences, Inc.

BMA 3F, 1-5-5 Minatojima-Minamimachi, Chuo-ku, Kobe 650-0047 JAPAN

TEL: 078-302-7091 / FAX: 078-302-7086

E-mail: info@carnabio.com

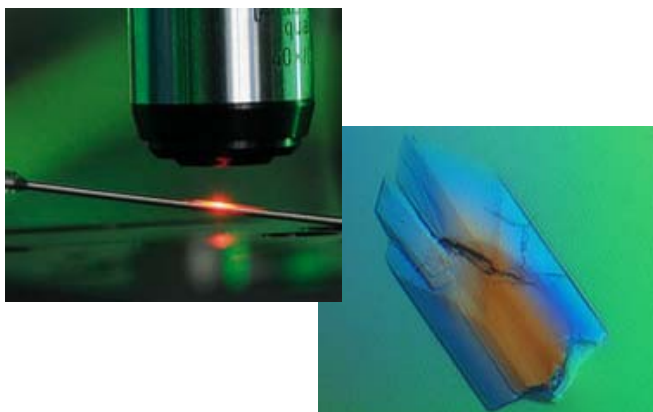
Co-Crystallography Service & Crystal Grade Proteins "XPRESS"

New Tools and Crystallography Services for Structure Based Drug Discovery

The Protein Kinases listed below are currently available for Co-Crystallography Services, and most(*1) can be offered as Crystal Grade Proteins(*2).

Available Targets - Protein Kinases

ABL	IRAK4
ACVR2A NEW	JAK1
ACVR2B NEW	JAK2
ALK	JAK3
ALK5	JNK1
ASK1	JNK2
AurA	JNK3
AXL1	LCK
AXL2 (MER)	LYN
B-RAF	MAPK11
CDK2	MET
C-FMS	MPS1
CHK1	NEK2
CK1 γ	p38
C-KIT	PAK1
DAPK3	PAK4
EGFR	PDK1
EPHB4	PGK1
ERK2	PIM1
ERK3/MAPK6	PIM2
FAK	PKA
FES	PKB
FGFR1 NEW	PKC ι NEW
FGFR2 NEW	PLK1
FGFR3 NEW	PLK4
FLT3	ROCK2
GSK3 β	RSK1
HASPIN	RSK2
HCK	SRC
IGF1-R	SYK



Courtesy of CRELUX GmbH

Available non-Kinase Targets

Chaperones	
DJ1	
HSP70	
HSP70/Bag1	
HSP90	
HSP90/cdc37	
Histone Acetyltransferase	
MYST1	
Histone Deacetylases	
HDAC4	
HDAC7	
HDAC8	
Histone Demethylases	
JMJD2A	
JMJD2C	
JMJD2D	
Histone Methyl-transferase	
EHMT1	
NHR	
ER α	
GR	
MCR	
PPAR α	
PPAR γ	
PPAR δ	
PR	
ROR γ	
Phosphatases	
PTP-1B	
PTPN22	
Protease	
Caspase3	
Poy (ADP-ribose) polymerases	
PARP1	
PARP3	
PARP12	
PARP14	
PARP15	
PARP5 (Tankyrase1)	
Sirtuin	
SIRT5	

(*1) Some proteins may require compounds at the time of expression, therefore, not available for individual purchase.

(*2) Proteins may not be active, crystallization is prioritized.

Visit our website and check the latest available targets!