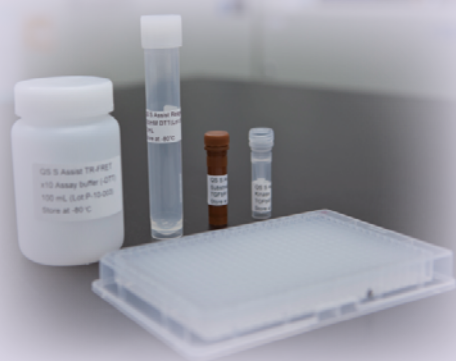


QuickScout Screening Assist™ Kits



QuickScout Screening Assist™ Kits are designed to speed you through compound screening, particularly secondary- and counter- screening operations, by providing necessary reagents and detailed assay protocols, for more than 300 human kinases.

After initial kit purchase, components may be purchased separately.

Minimum order requirements apply.

* The dilution ratio and MSA measuring parameters are different by kinase.

Advantages of Carna's Assay Kits...

- Prepared utilizing the extensive expertise of our profiling services
- Ready-To-Run products & protocols
- One kit allows you multiple assays

Designed for primary, in-house screening procedures - Lead Generation through Lead Optimization!

* As of June 21, 2018

Platform	Minimum Size	Available Target	Kit Components
Mobility Shift Assay QSS Assist™ MSA This MSA kit works best using LabChip® technology from PerkinElmer, Inc.	400dp Equivalent to 1 x 384-well plate	295 Kinases	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Metal included) ● Assay Buffer ● Termination Buffer ● Assay Protocol (Separation conditions included) * Protocol sample is available online.
FP (IMAP™) QSS Assist™ FP	800dp Equivalent to 2 x 384-well plate	81 Kinases	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Assay Protocol
TR-FRET QSS Assist™ TR-FRET	800dp Equivalent to 2 x 384-well plate	28 Kinases	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Assay Protocol
ELISA QSS Assist™ ELISA	100dp Equivalent to 1 x 96-well plate	32 Kinases	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Antibody for ELISA (except for TTK & WEE1) ● Assay Protocol
ADP-Glo™ QSS Assist™ ADP-Glo™	400dp 1 x 384-well plate 相当	11 Kinases	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Solution ● Kinase dilution Buffer (Metal included) ● Assay Buffer ● MgCl2 solution for detection reagent ● Assay Protocol

Each kinase kit is made-to-order, with turnaround time of 2-3 weeks.

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

Carna's wholly-owned subsidiary

CarnaBio USA, Inc.

209 West Central Street, Suite 307, Natick, MA 01760 USA

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

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