

# 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. Each assay is designed to be completed in 2-3 hours at ATP levels approximating Km value. Each kinase kit is made-to-order, with turnaround time of 2-3 weeks.

## 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
- Up to 4 kinases can be selected per kit (MSA 400dp only)



**Time & Cost Effective!**

## READY-TO-RUN Products & Protocols for 312 Kinase Assays

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

### Mobility Shift Assay (MSA) QSS Assist™ MSA

This MSA kit works best using LabChip® technology from Caliper Life Sciences, Inc.



**800dp Set**  
(Equivalent to 2 x 384-well plate)

**272 Kinases**

ONE kinase may be selected

**400dp Set**  
(Equivalent to 1 x 384-well plate)

**209 Kinases**

Select either 1, 2, or 4 kinases

### QSS Assist™ MSA Kit Components

- Protein Kinase
- Substrate Mixture (ATP, Metal included)
- Assay Buffer
- Termination Buffer
- Assay Protocol (Separation conditions included) ← Protocol sample is available online.

After initial kit purchase, components may be purchased separately. Minimum order requirements apply. For more details, please visit our website.

### FP (IMAP™) QSS Assist™ FP

**77 Kinases**

**Kit Components**

- Protein Kinase
- Substrate Mixture
- Assay Buffer
- Assay Protocol

800dp/set (Equivalent to 2 x 384-well plate)

### TR-FRET QSS Assist™ TR-FRET

**22 Kinases**

**Kit Components**

- Protein Kinase
- Substrate Mixture
- Assay Buffer
- Assay Protocol

800dp/set (Equivalent to 2 x 384-well plate)

### ELISA QSS Assist™ ELISA

**110 Kinases**

**Kit Components**

- Protein Kinase
- Substrate Mixture
- Assay Buffer
- Assay Protocol

500dp/set (Equivalent to 5 x 96-well plate)

September 25, 2012 Updated

## 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

