



## ADP-Glo™ Kinase Profiling Application Notes

# CMGC KINASE KSPS: CMGC-1

# Kinase Selectivity Profiling System: CMGC-1

By Jacquelyn Hennek, M.S., Said A. Goueli, Ph.D., and Hicham Zegzouti, Ph.D., Promega Corporation

## Scientific Background:

Kinase Selectivity Profiling System CMGC-1 is a set of kinases from the CMGC Kinase Family presented in an easy to use 8-tube strip format. When diluted, the kinase stock volumes are standardized to generate optimal ATP to ADP conversion with a signal to background ratio over 10-fold when their activities are detected using the ADP-Glo™ Kinase Assay (Fig. 1). The substrate stocks are standardized in a similar fashion and are located in a second strip at corresponding positions. Kinase Selectivity Profiling Systems can be used to generate single-dose inhibitor selectivity profiles for as many inhibitors as desired (Fig. 2A) or to study dose response curves for an inhibitor (Fig. 2B).

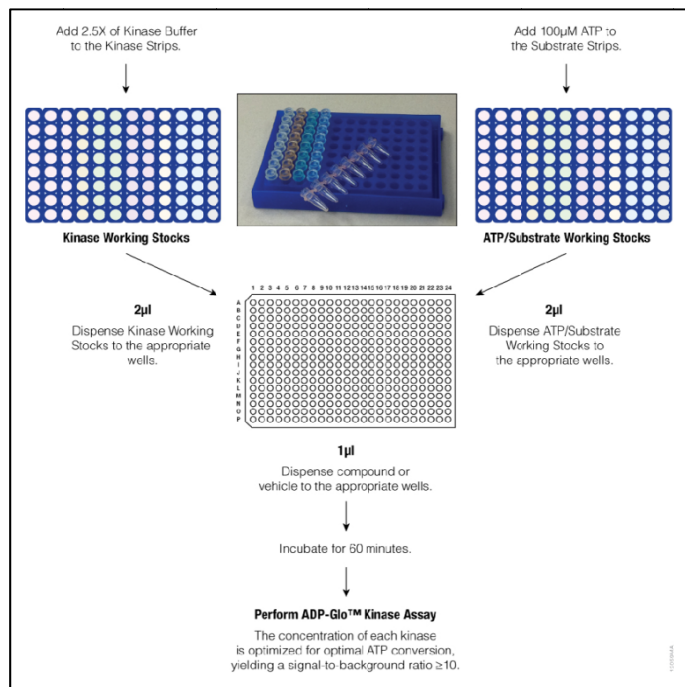
CMGC-1		
CMGC Family		
	Kinase Strip	Substrate Strip
A	ERK2	MBP
B	GSK3β	GSK3 Substrate
C	JNK1	p38 Substrate
D	JNK3	p38 Substrate
E	p38α	p38 Substrate
F	p38β	p38 Substrate
G	p38δ	p38 Substrate
H	p38γ	p38 Substrate

## ADP-Glo™ Kinase Assay

### Description

ADP-Glo™ Kinase Assay is a luminescent kinase assay that measures ADP formed from a kinase reaction; ADP is converted into ATP, which is converted into light by Ultra-Glo™ Luciferase.

The luminescent signal positively correlates with ADP amount and kinase activity. The assay is well suited for measuring the effects chemical compounds have on the activity of a broad range of purified kinases—making it ideal for both primary screening as well as kinase selectivity profiling (Fig. 2).



**Figure 1. Kinase Selectivity Profiling System Overview.** Kinases are provided at either 25X or 50X concentrations in an 8-tube strip, and substrates/cofactors are provided at 3.3X concentrations in a separate 8-tube strip. One-step dilutions directly into the strips produce sufficient Kinase and ATP/Substrate Working Stocks for 25 kinase reactions. Kinase reactions are performed using 1µL of compound, 2µL of Kinase Working Stock, and 2µL of ATP/Substrate Working Stock. After 1 hour incubation at room temperature, kinase activity is quantified using the ADP-Glo™ Kinase Assay. The luminescent signal generated by the ADP-Glo™ Kinase Assay is proportional to ADP concentration and correlated with kinase activity.

For detailed protocols on strip preparation, single-dose inhibition profiles, and creating dose-response curves, see *The Kinase Selectivity Profiling System* Technical Manual #TM421, available at [www.promega.com/protocols/tm421](http://www.promega.com/protocols/tm421)



### Preparation of Kinase and ATP/Substrate Working Stocks:

- Add 95µl of 2.5X Kinase Buffer to all tubes in the Kinase Strip.
- Add 15µl of 100µM ATP to all tubes in the Substrate/Cofactors Strip.

### Single-Dose Inhibition Profile:

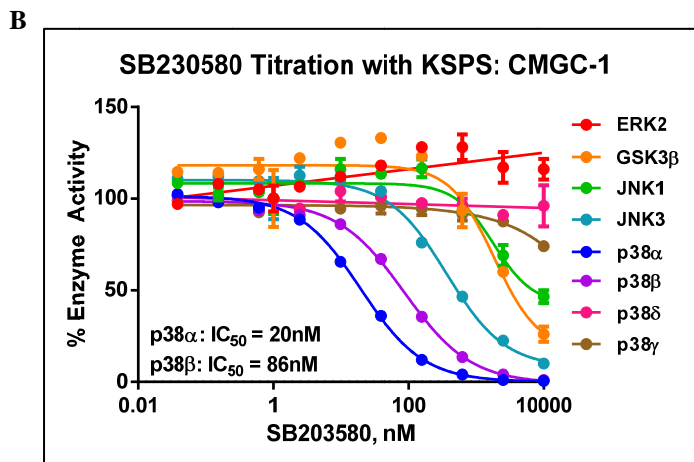
- Setup Kinase Reactions and No Compound Controls:
  - 1µl of compound or vehicle (5% DMSO)
  - 2µl of Kinase Working Stock
- Setup No Kinase Controls:
  - 1µl vehicle (5% DMSO)
  - 2µl of Kinase Buffer
- Incubate at room temperature for 10 minutes.
- Add 2µl of ATP/Substrate Working Stock.
- Incubate at room temperature for 60 minutes.
- Perform ADP detection using ADP-Glo™ Kinase Assay.

### Dose-Response Curves:

- Setup Kinase Reactions:
  - 1µl of 5X compound serial dilution
  - 2µl of Kinase Working Stock
- Setup No Kinase Controls:
  - 2µl of Kinase Buffer in place of Kinase Working Stock
- Incubate at room temperature for 10 minutes.
- Add 2µl of ATP/Substrate Working Stock.
- Incubate at room temperature for 60 minutes.
- Perform ADP detection using ADP-Glo™ Kinase Assay.

**A**

		SB203580	Roscovitine
<b>KSPS: CMGC-1</b>	<b>ERK2</b>	97	99
	<b>GSK3β</b>	55	96
	<b>JNK1</b>	84	107
	<b>JNK3</b>	30	106
	<b>p38α</b>	3	101
	<b>p38β</b>	10	101
	<b>p38δ</b>	105	107
	<b>p38γ</b>	94	104



**Figure 2. KSPS: CMGC-1 Profiling Data.** (A) KSPS: CMGC-1 kinase activities were determined in the presence of 1µM SB203580 or Roscovitine. % Activity values were calculated using No Compound and No Kinase Controls and are shown above. Red < 20%; White 20-60%; Blue > 60%. (B) SB203580 dose response curves were created with KSPS: CMGC-1 to determine the potency (IC<sub>50</sub>) and selectivity of the inhibitor. IC<sub>50</sub> values are comparable to literature values<sup>(1)</sup>.

<sup>(1)</sup> Cuenda, A. et. al.; FEBS Letters. 1995, 364, 229.

### Assay Components and Ordering Information:



#### Products

ADP-Glo™ Kinase Assay  
 Kinase Selectivity Profiling System: CMGC-1  
 Kinase Selectivity Profiling System: CMGC-1 + ADP-Glo™ Assay

#### Company

Promega  
 Promega  
 Promega

#### Cat.#

V6930  
 V6854  
 V6855

Kinase Buffer: 40mM Tris, pH 7.5; 20mM MgCl<sub>2</sub>; 0.1mg/ml BSA; 50µM DTT.