

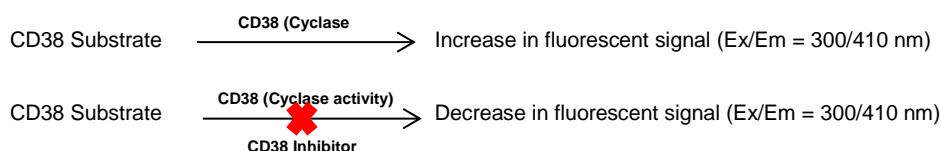
## CD38 (Cyclase) Inhibitor Screening Kit (Fluorometric)

09/21

(Catalog # K2102-100; 100 assays; Store at -20 °C)

### I. Introduction:

Cluster of differentiation 38 (CD38), also known as cyclic ADP ribose hydrolase or ADP-ribosyl cyclase is a multifunctional enzyme that catalyzes the synthesis and hydrolysis of Nicotinamide Adenine Dinucleotide (NAD) to Nicotinamide and ADP-ribose (ADPR). Additionally, CD38 generates second messengers, cyclic ADP-ribose using NAD as the substrate. Due to its role in NAD<sup>+</sup> metabolism, the study of CD38 and its functions has been of great importance. It is found on the surface of many immune cells, including plasma B cells, natural killer cells, CD4<sup>+</sup>, CD8<sup>+</sup>, etc. Elevated levels of CD38 are associated with aging, obesity, diabetes, heart disease, asthma, inflammation and tumorigenesis etc. **BioVision's CD38 (Cyclase) Inhibitor Screening Kit** is a plate-based fluorometric assay designed to screen, study and characterize potential inhibitors of CD38 cyclase activity. The assay utilizes a selective CD38 substrate to generate a fluorescent signal measured at Ex/Em = 300/410 nm. In the presence of potential CD38 cyclase inhibitors, the fluorescent signal is reduced. The assay is quick, easy, and sensitive for high-throughput screening of CD38 inhibitors. Additionally, the kit includes a CD38 Inhibitor as a control inhibitor.



### II. Application:

- Screening or characterizing CD38 (Cyclase) inhibitors.

### III. Kit Contents:

Components	K2102-100	Cap Code	Part Number
CD38 Assay Buffer	25 ml	WM	K2102-100-1
CD38 Substrate	1 vial	White	K2102-100-2
CD38, Human Recombinant	1 vial	Blue	K2102-100-3
CD38 Inhibitor	50 $\mu$ l	Yellow	K2102-100-4

### IV. User Supplied Reagents and Equipment:

- 96-well white plate with flat bottom
- Multi-well spectrophotometer

### V. Storage Conditions and Reagent Preparation:

Store kit at -20 °C, protected from light. Briefly centrifuge all small vials prior to opening. Read the entire protocol before performing the assay.

- **CD38 Assay Buffer:** Store at 4 °C or -20 °C. Bring to room temperature (RT) before use.
- **CD38 Substrate:** Reconstitute the vial in 220  $\mu$ l dH<sub>2</sub>O. Divide into aliquots and store at -20 °C. Keep on ice during use.
- **CD38, Human Recombinant:** Reconstitute the vial in 110  $\mu$ l CD38 Assay Buffer. Divide into aliquots and store at -20 °C. Keep on ice during use. Avoid repeated freeze-thaw cycles.
- **CD38 Inhibitor (5 mM in DMSO):** Warm to RT. Divide into aliquots and store at -20 °C.

### VI. CD38 Inhibitor Screening Protocol:

**1. CD38, Human Recombinant Enzyme Dilution:** Prepare 1:5 dilution of the CD38 enzyme using CD38 Assay Buffer. Mix thoroughly and keep on ice. Add 2.5  $\mu$ l of diluted CD38 enzyme into the desired wells of a 96-well white plate labeled as **Sample**, **Solvent Control**, **Inhibitor Control** and **Enzyme Control**. Adjust the volume of all wells to 25  $\mu$ l using CD38 Assay Buffer.

**2. Screening Test Inhibitor(s):** Dissolve the Test Inhibitor(s) in an appropriate solvent to make 100X stock solution. Dilute the stock Test Inhibitor to 4X using CD38 Assay Buffer. Add 25  $\mu$ l of diluted Test Inhibitor into the **Sample** well(s). Add 25  $\mu$ l of 4X Solvent (4X final well solvent concentration) into the **Solvent Control** well. **Note:** Solvents used to solubilize the Test Inhibitor(s) might affect the enzymatic activity. Thus, prepare a **Solvent Control** well by adding 25  $\mu$ l of solution with the same final concentration of solvent in assay buffer that is used to dissolve the Test Inhibitor(s).

**3. Enzyme Control, Background Control and Inhibitor Control Preparation:** Add 25  $\mu$ l of CD38 Assay Buffer to the **Enzyme Control** well. For **Background Control**, add 50  $\mu$ l of CD38 Assay Buffer in a separate well. To the **Inhibitor Control** well, add 2  $\mu$ l of 5 mM CD38 Inhibitor and adjust the volume to 50  $\mu$ l/well by adding 23  $\mu$ l CD38 Assay Buffer. At this stage, the volume of all wells including Sample, Solvent Control, Inhibitor Control, Enzyme Control and Background Control is 50  $\mu$ l/well.

**IC<sub>50</sub> estimation (Optional):** Prepare several dilutions of the Test Inhibitor(s) in CD38 Assay Buffer while maintaining the consistent final Solvent Concentration in all wells. Add 25  $\mu$ l of each dilution into the designated wells.

**4. CD38 Substrate Mix Preparation:** Mix enough CD38 Substrate Mix for the number of assays to be performed. For each well, prepare 50  $\mu$ l CD38 Substrate Mix containing

	<u>CD38 Substrate Mix</u>
CD38 Assay Buffer	48 $\mu$ l
CD38 Substrate	2 $\mu$ l

Add 50  $\mu$ l CD38 Substrate Mix to Sample, Solvent Control, Inhibitor Control, Enzyme Control and Background Control wells and mix well. The total reaction volume is 100  $\mu$ l/well.

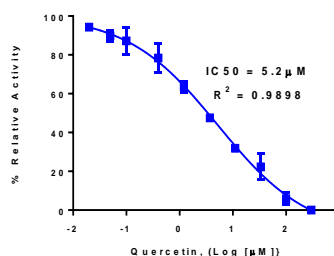
**5. Measurement:** Measure the fluorescence in kinetic mode (Ex/Em = 300/410 nm) at 37 °C for 30-60 min. Choose any two time points ( $t_1$  &  $t_2$ ) in the linear range of the plot and obtain the corresponding RFU values.

**Note:** The Enzyme progressive curve is hyperbolic, with an initial linear portion followed by progressively slower reaction. Use the initial portion to check the linear range of the reaction.

**6. Calculation:** Subtract the RFU of the BC well from all Test Inhibitor(s) [S], Enzyme Control [EC], Solvent Control [SC] and Inhibitor Control [IC] wells. Obtain  $\Delta$ RFU for S, EC, SC and IC by subtracting RFU at time  $t_1$  from RFU at time  $t_2$ , such that  $t_2$  and  $t_1$  is within a linear range of the assay. If  $\Delta$ RFU of Solvent Control [SC] is significantly different from  $\Delta$ RFU of Enzyme Control [EC], use its values to determine the effect of test inhibitors dissolved in the same solvent.

Calculate the relative % inhibition of the Test Inhibitor(s) as below:

$$\% \text{ Relative Inhibition} = \frac{\Delta\text{RFU}[\text{EC}] - \Delta\text{RFU}[\text{S}]}{\Delta\text{RFU}[\text{EC}]} \times 100$$



**Figure:** Inhibition of CD38 cyclase activity by Quercetin.  $IC_{50}$  was calculated to be  $5.2 \pm 0.8 \mu\text{M}$ . Assay was performed following the kit protocol.

#### VII. Related Products:

Human CellExp™ CD38, human recombinant (Cat. # P1014-10, 50)  
 Human CellExp™ CD38, Mouse Recombinant (Cat. # P1338-10, 50)  
 CD38 (Cyclase) Activity Assay Kit (Fluorometric) (Cat. # K2042-100)  
 CD38 (Hydrolase) Activity Assay Kit (Fluorometric) (Cat. # K2095-100)  
 CD38 (Hydrolase) Inhibitor Screening Kit (Fluorometric) (Cat. # K2086-100)  
 NAD+/NADH Quantification Colorimetric Kit (Cat. # K337-100)  
 PicoProbe™ NADH Fluorometric Assay Kit (Cat. # K338-100)  
 EZScreen™ NAD+/NADH Colorimetric Assay Kit (384-well) (Cat. # K958-400)  
 Apigenin (Cat. # 2508-25, 100)  
 Quercetin, Dihydrate (Cat. # 1773-250, 1000)

**FOR RESEARCH USE ONLY! Not to be used on humans.**