

ab308152 – CD38 (Cyclase) Inhibitor Screening Kit (Fluorometric)

A 96-well plate based colorimetric assay to measure the activation of human c-Fos in nuclear extracts or cell lysates.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

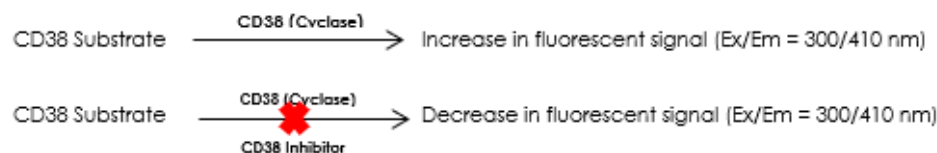
<http://www.abcam.com/ab308152>

Storage and Stability

The entire Assay kit may be stored at -20 °C and protected from light.

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⁺ 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⁺, CD8⁺, etc. Elevated levels of CD38 are associated with aging, obesity, diabetes, heart disease, asthma, inflammation and tumorigenesis etc. **The 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.



Materials Supplied

Item	Quantity
CD38 Assay Buffer	25 mL
CD38 Substrate	1 vial
CD38, Human Recombinant	1 vial
CD38 Inhibitor	50 μ l

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

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

Reagent Preparation

- Store the 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 μ l dH₂O. Divide into aliquots and store at -20 °C. Keep on ice during use.

CD38, Human Recombinant: Reconstitute the vial in 110 μ 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.

Sample Preparation

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 μ 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 μ l using CD38 Assay Buffer.

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 μ l of diluted Test Inhibitor into the **Sample** well(s). Add 25 μ 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 μ l of solution with the same final concentration of solvent in assay buffer that is used to dissolve the Test Inhibitor(s).

Enzyme Control, Background Control and Inhibitor Control Preparation: Add 25 μ l of CD38 Assay Buffer to the Enzyme Control well. For Background Control, add 50 μ l of CD38 Assay Buffer in a separate well. To the Inhibitor Control well, add 2 μ l of 5 mM CD38 Inhibitor and adjust the volume to 50 μ l/well by adding 23 μ 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 μ l/well.

IC50 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 μ l of each dilution into the designated wells.

Assay Procedure

1. **CD38 Substrate Mix Preparation:** Mix enough CD38 Substrate Mix for the number of assays to be performed. For each well, prepare 50 µl CD38 Substrate Mix containing

	<u>CD38 Substrate Mix</u>
CD38 Assay Buffer	48 µl
CD38 Substrate	2 µl

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

2. **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.
3. **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 Δ 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 Δ RFU of Solvent Control [SC] is significantly different from Δ 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:

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

For technical support contact information, visit: www.abcam.com/contactus