



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### **CD38 Inhibitor Screening Assay Kit (Cyclase Activity)**

**Catalog # 71275**

**BACKGROUND :** CD38, a differentiation antigen of B lymphocytes, is a type II integral membrane protein. It is also known as ADP-ribosyl cyclase and nicotinamide adenine dinucleotide (NAD<sup>+</sup>) glycohydrolase. Through its production of cyclic ADP-ribose, CD38 modulates calcium-mediated signal transduction in various cells, including pancreatic  $\beta$  cells. CD38 is a prognostic biomarker for acute B lymphoblastic leukemia.

**DESCRIPTION:** The *CD38 Inhibitor Screening Assay Kit (Cyclase Activity)* is designed to measure the cyclase activity of CD38 for screening and profiling applications. The CD38 assay kit comes in a convenient 96-well format, with purified recombinant CD38 enzyme, its substrate nicotinamide guanine dinucleotide (NGD<sup>+</sup>), and CD38 assay buffer for 100 enzyme reactions. In addition, the kit includes the CD38 inhibitor quercetin for use as a positive control.

#### **COMPONENTS:**

Catalog #	Reagent	Amount	Storage	
71277	CD38, His-Tag (Human), HiP™	25 $\mu$ g	-80°C	<b>Avoid multiple freeze/thaw cycles!</b>
	3x CD38 assay buffer	4 ml	-20°C	
	CD38 substrate NGD <sup>+</sup>	50 $\mu$ l	-20°C	
	Quercetin (50 mM DMSO)	100 $\mu$ l	-20°C	
79685	Black 96-well plate	1	Room Temp.	

#### **MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Adjustable micropipettor and sterile tips  
Fluorescent microplate reader  
Rotating or rocker platform

**APPLICATIONS:** Great for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

**STABILITY:** Up to 6 months from date of receipt, when stored as recommended.

**REFERENCE:** Wei, W., *et al.*, *World J. Biol. Chem.* 2014 **5**(1):58-67.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

1. Thaw **3x CD38 assay buffer** on ice.
2. Prepare the master mixture (20  $\mu$ l per well): N wells x (10  $\mu$ l **3x CD38 assay buffer** + 10  $\mu$ l water). Add 20  $\mu$ l to every well.

	Positive Control	Test Inhibitor	Blank
3x CD38 assay buffer	10 $\mu$ l	10 $\mu$ l	10 $\mu$ l
Water	10 $\mu$ l	10 $\mu$ l	10 $\mu$ l
Test Inhibitor	-	10 $\mu$ l	-
Inhibitor Buffer (no inhibitor)	10 $\mu$ l	-	10 $\mu$ l
1x CD38 assay buffer	-	-	15 $\mu$ l
CD38 (16.7 ng/ $\mu$ l)	15 $\mu$ l	15 $\mu$ l	-
NGD <sup>+</sup> (diluted)	5 $\mu$ l	5 $\mu$ l	5 $\mu$ l
Total	50 $\mu$ l	50 $\mu$ l	50 $\mu$ l

3. Add 10  $\mu$ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the wells labeled "Positive Control" and "Blank", add 10  $\mu$ l of the same solution without inhibitor (Inhibitor buffer).
4. Prepare **1x CD38 assay buffer** by diluting **3x CD38 assay buffer** with water. Dilute only enough buffer required for the assay. Store remaining **3x CD38 assay buffer** at -20°C in single-use aliquots. For 100 reactions, prepare 6 ml **1x CD38 assay buffer** by mixing 2 ml of **3x CD38 assay buffer** with 4 ml water.
5. To the wells designated as "Blank", add 15  $\mu$ l of **1x CD38 assay buffer**.
6. Thaw **CD38** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Calculate the amount of **CD38** required for the assay and dilute enzyme to 16.7 ng/ $\mu$ l with **1x CD38 assay buffer** (250 ng/well). Aliquot remaining **CD38** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: CD38 enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
7. Add 15  $\mu$ l of diluted **CD38** enzyme to the wells designated "Positive Control" and "Test Inhibitor Control". Cover the plate and incubate 1 hour at room temperature with slow shaking.

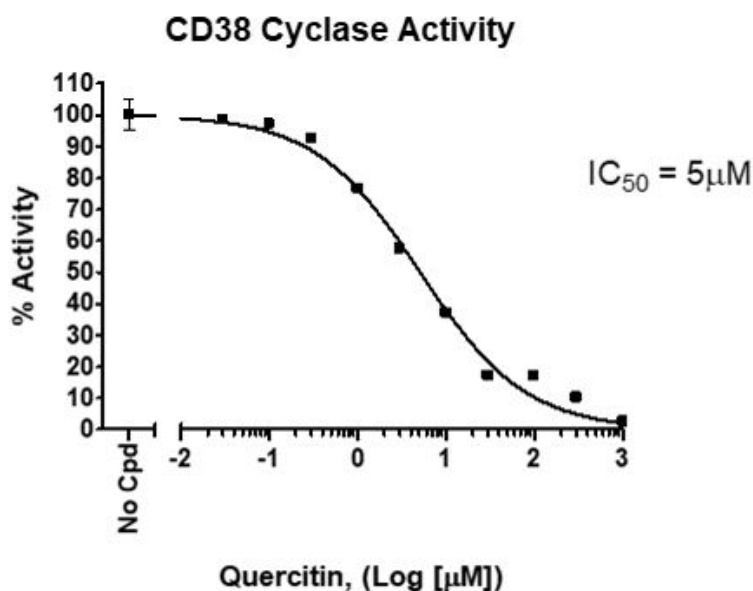
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

8. During incubation, dilute **NGD<sup>+</sup>** 10-fold with **1x CD38 assay buffer**. Dilute only the amount required for the assay. Store remaining **NGD<sup>+</sup>** at -20°C in single use aliquots. Discard any unused diluted **NGD<sup>+</sup>** after use.
9. After the 1 hour incubation, remove the plate and add 5 µl of diluted **NGD<sup>+</sup>**.
10. Place plate into plate-reading fluorimeter and prepare to measure.
11. After 4-6 minutes, measure the plate using a fluorimeter capable of excitation at 300 nm and detection of emitted light at 410 nm. The "Blank" value is subtracted from all other values.

**Example of Assay Results:**

CD38 inhibition by quercetin, measured using the **CD38 Inhibitor Screening Assay Kit (cyclase activity)**, BPS Bioscience Cat. # 71275. Fluorescence was measured using a Bio-Tek microplate reader. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<b><u>Product Name</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
CD38, His-Tag (Human), HiP™	71227	100 µg
5'-Nucleotidase/CD73, His-tag	71184	50 µg
Quercetin	27214	5 g
CD73 Inhibitor Screening Assay Kit	72055	96 rxns
CD73 Inhibitor Screening Assay Kit	72058	384 rxns
Adenosine Deaminase (ADA), His-tag	70016	100 µg
NAD+, Biotin-Labeled	80610	500 µl
NAMPT (PBEF1)	71098	50 µg
NAMPT (PBEF1)	91004	50 µg
NMNAT, His-tag	71090	100 µg
TCF/LEF Reporter Kit	60500	500 rxns
TCF/LEF reporter-HEK293 cell line	60501	2 vials

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)