



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data sheet
CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)
Catalog #79672
Size: 384 reactions

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) glycohydrolase. Through its production of cyclic ADP-ribose, CD38 modulates calcium-mediated signal transduction in various cells, including pancreatic β cells. The major enzymatic activity of CD38 is the hydrolysis of NAD. CD38 is a prognostic biomarker for acute B lymphoblastic leukemia.

DESCRIPTION: The *CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)* is designed to measure the glycohydrolase activity of CD38 for screening and profiling applications. The CD38 assay kit comes in a convenient high-throughput 384-well format, with purified recombinant CD38 enzyme, its substrate N6-etheno-NAD (ϵ -NAD), and CD38 assay buffer for 384 enzyme reactions. In addition, the kit includes the CD38 inhibitor apigenin for use as a control inhibitor.

COMPONENTS:

Catalog #	Component	Amount	Storage	
71277	CD38, His-Tag (Human), HiP™	12 μ g	-80 °C	Avoid multiple freeze/thaw cycles!
	4x CD38 hydrolase buffer	10 ml	-20 °C	
	CD38 substrate (ϵ -NAD)	200 μ l	-20 °C	
	Apigenin (50 mM DMSO)	10 μ l	-20 °C	
79961	Black 384-well plate			

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.

REFERENCES:

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

1. Thaw **4x CD38 hydrolase buffer** on ice.
2. Prepare the master mixture (10 μ l per well): N wells x (5 μ l **4x CD38 hydrolase buffer** + 5 μ l water). Add 10 μ l to every well.

	Positive Control	Negative Control	Test Inhibitor	Blank
4x CD38 hydrolase buffer	5 μ l	5 μ l	5 μ l	5 μ l
Water	5 μ l	5 μ l	5 μ l	5 μ l
Test Inhibitor	-	-	10 μ l	
Inhibitor buffer (no inhibitor)	10 μ l	-	-	10 μ l
Apigenin	-	10 μ l	-	-
1x CD38 hydrolase buffer	-	-	-	20 μ l
CD38 (1.5 ng/ μ l)	20 μ l	20 μ l	20 μ l	-
ϵ -NAD (diluted)	10 μ l	10 μ l	10 μ l	10 μ l
Total	50 μl	50 μl	50 μl	50 μl

3. Prepare **1x CD38 hydrolase buffer** by diluting **4x CD38 hydrolase buffer** with water. Dilute only enough buffer required for the assay. Store remaining **4x CD38 hydrolase buffer** at -20°C in single-use aliquots. For 384 reactions, prepare 10 ml **1x CD38 hydrolase buffer** by mixing 2.5 ml of **4x CD38 hydrolase buffer** with 7.5 ml water.
4. Add 10 μ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the wells labeled "Positive Control" and "Blank", add 10 μ l of the same solution without inhibitor (Inhibitor buffer).
*Optional: For the well labeled "Negative Control", add 10 μ l apigenin, diluted 0.1 – 100 μ M in 1x CD38 hydrolase buffer.
5. To the wells designated as "Blank", add 20 μ l of **1x CD38 hydrolase 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 1.5 ng/ μ l with **1x CD38 assay buffer** (30 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.

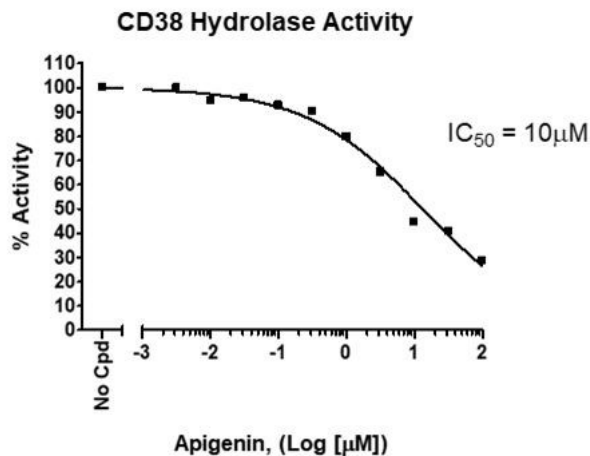
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Please visit our website at: www.bpsbioscience.com

7. Add 20 μ l of diluted **CD38** enzyme to the wells designated "Positive Control", "Negative Control", and "Test Inhibitor Control". Cover the plate and incubate 30 minutes at room temperature with slow shaking.
8. During incubation, dilute ϵ -**NAD** 20-fold with **1x CD38 hydrolase buffer**. Dilute only the amount required for the assay. Store remaining ϵ -**NAD** at -20°C in single use aliquots. Discard any unused diluted ϵ -**NAD** after use.
9. After the 30 minutes incubation, remove the plate and add 10 μ l of diluted ϵ -**NAD**.
10. Place plate into plate-reading fluorimeter and prepare to measure.
11. After 10 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 apigenin, measured using the **CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)**, BPS Bioscience #79672. 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

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog#</u>	<u>Size</u>
CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79287	96 rxns
CD38, His-Tag (Human), HiP™	71227	100 µg
CD39, His-tag	71284	20 µg
CD73, Avi, His-tag (Mouse)	72523	100 µg
CD73, His-tag	71184	50 µg
CD73 Inhibitor Screening Assay Kit	72055	96 rxns
CD73 Inhibitor Screening Assay Kit	72058	384 rxns
CD38 Inhibitor Screening Assay Kit	71275	96 rxns
CD39 Inhibitor Screening Assay Kit	79278	96 rxns
Quercetin	27214	5 g
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

Please visit our website at: www.bpsbioscience.com