



6405 Mira Mesa Blvd Ste. 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Data Sheet

IDH1(R132H) Assay Kit

Catalog # 79376

BACKGROUND : Isocitrate dehydrogenases are enzymes that catalyze the oxidative decarboxylation of isocitrate to 2-oxoglutarate. Mutations in isocitrate dehydrogenases 1 and 2 (IDH1 and IDH2) have been reported in a multitude of human cancers. While the wild type IDH1, produces α -ketoglutarate (α -KG) through the reduction of NADP⁺ to NADPH, the IDH1(R132H) mutant catalyzes conversion of α -ketoglutarate to 2-hydroxyglutarate (2-HG) by oxidizing NADPH to NADP⁺. Genomic studies have identified the IDH1(R132H) mutation in various cancer types, including glioma, chondrosarcoma, and AML as well as a small percentage of prostate, lung and colon cancers.

DESCRIPTION: The *IDH1(R132H) Assay Kit* is designed to measure IDH1(R132H) activity for screening and profiling applications by measuring NADPH consumption. The *IDH1(R132H) Assay Kit* comes in a convenient 96-well format, with enough purified recombinant IDH1(R132H) enzyme, α -KG, NADPH, NADPH detection reagents, and assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	Avoid multiple freeze/thaw cycles!
71099	IDH1(R132H)	5 μ g	-80°C	
	1x IDH1 Assay Buffer	20 ml	-20°C	
	α -Ketoglutarate (40 mM)	1 ml	-20°C	
	NADPH (500 μ M)	2 x 100 μ l	-20°C	
	NADPH Detection Reagent A	2 x 1 mg	-20°C	
	NADPH Detection Reagent B	2 x 1 mg	-20°C	
79685	96-well plate, black	2	Room Temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

NADPH (optional)*

Microplate reader capable of reading Fluorescence

Adjustable micropipettor and sterile tips

**Note: NADPH is not stable for multiple freeze/thaw cycles. We have provided the reagent as 2 x 100 μ l to allow you to use only half the plate at a time. If you plan to use the same plate for additional experiments, please purchase additional NADPH and prepare a fresh 500 μ M stock in H₂O at the time of the assay.*

APPLICATIONS: Useful for screening small molecular inhibitors for drug discovery and HTS applications.

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: support@bpsbioscience.com



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

STABILITY: Up to 6 months when stored as recommended but varies by components.

REFERENCE: Wang, F., *et. al. Science* **340**:622-626 (2013)

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Thaw **1x IDH1 Assay Buffer**, **α -Ketoglutarate (40 mM)**, and **NADPH (500 μ M)**.
- 2) Prepare the master mixture (70 μ l per well): N wells x (58 μ l **1x IDH1 Assay Buffer**, + 2 μ l **NADPH (500 μ M)**, and + 10 μ l **α -Ketoglutarate (40 mM)**). Add 70 μ l to every well.

	Positive Control	Test Inhibitor	Blank
1x IDH1 Assay buffer	58 μ l	58 μ l	58 μ l
NADPH (500 μ M)	2 μ l	2 μ l	2 μ l
α -Ketoglutarate (40 mM)	10 μ l	10 μ l	10 μ l
Test Inhibitor	–	10 μ l	–
Inhibitor Buffer (no inhibitor)	10 μ l	–	10 μ l
1x IDH1 Assay buffer	–	–	20 μ l
IDH1(R132H) enzyme; 2-2.5 ng/ μ l~2.5 ng/ μ l	20 μ l	20 μ l	–
Total	100 μ l	100 μ l	100 μ l

- 3) Add 10 μ l of Inhibitor solution of each well labeled as "Test Inhibitor." For the "Positive Control" and "Blank," add 10 μ l of the same solution without inhibitor (Inhibitor buffer).

Note: Some inhibitors require preincubation with IDH1(R132H) to exhibit maximum effect. Also, depending on the mechanism of the inhibitor, the pre-incubation may require adding one or both substrates (α -Ketoglutarate and NADPH) separately (step 2). For example, for IDH-C100 inhibition, we preincubate with IDH1(R132H) enzyme for 90 min in the presence of NADPH, and then initiate the reaction by adding the α -Ketoglutarate.

- 4) To the wells designated as "Blank," add 20 μ l **1x IDH1 Assay buffer**.

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: support@bpsbioscience.com



6405 Mira Mesa Blvd Ste. 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

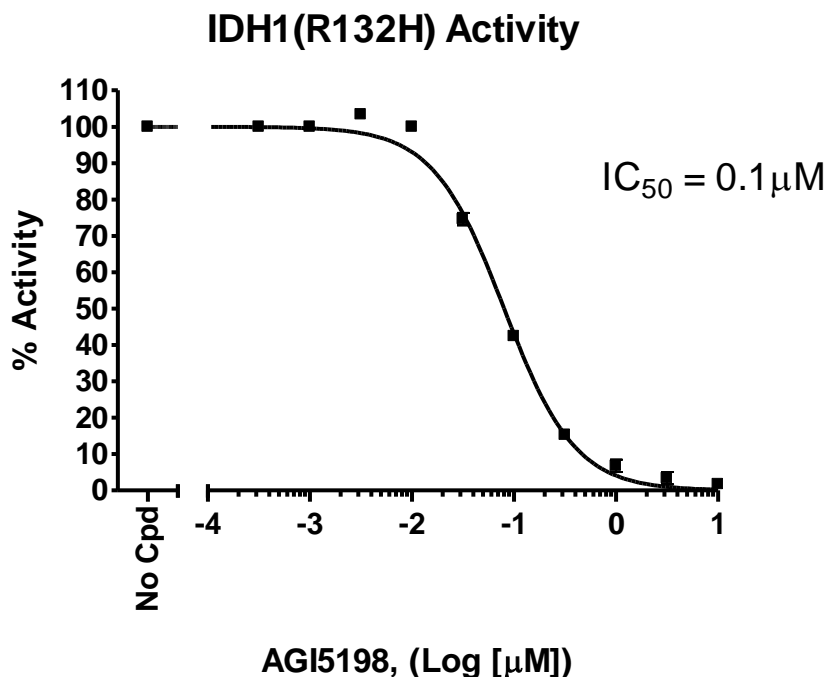
- 5) Thaw **IDH1(R132H)** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of **IDH1(R132H)** required for the assay and dilute enzyme to 2~2.5 ng/μl for **IDH1(R132H)** with **1x IDH1 Assay Buffer**. Store remaining undiluted enzyme in aliquots at -80°C. *Note: IDH1(R132H) enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 6) Initiate reaction by adding 20 μl of diluted **IDH1(R132H)** enzyme to the wells designated "Positive Control" and "Test Inhibitor." Incubate at room temperature for 1 hour.
- 7) Immediately before terminating the reaction, add 1 ml distilled water to **NADPH Detection Reagent A** and 0.4 ml water to **NADPH Detection Reagent B** and mix well to completely dissolve the reagents. Store the tubes on ice.
- 8) Prepare the NADPH detection reagent by adding 180 μl of **NADPH Detection Reagent A** and 10 μl of **NADPH Detection Reagent B** to 4,810 μl of **1x IDH1 Assay Buffer**. *Note: NADPH detection reagents are not stable so it must be used immediately; discard any unused reagent. The kit provides enough reagents for two separate assays.*
- 9) Add 25 μl of NADPH detection reagent to the wells of a new plate (provided). Using a multipipettor, remove 50 μl of the **IDH1(R132H)** reaction mixture and add to the wells containing the NADPH detection reagent. Cover the plate with aluminum foil and incubate for 5 minutes at room temperature.
- 10) Read the fluorescence at Ex544 nm/Em600 nm using a microplate reader.

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: support@bpsbioscience.com

Example of Assay Results:



Inhibition of IDH1(R132H) by AGI-5198, measured using the *IDH1(R132H) assay kit* (Cat. #79376). For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
IDH2, FLAG-tag	71074-1	50 μg
IDH2, FLAG-tag	71074-2	100 μg
IDH2 (R140Q), FLAG-tag	71100-1	50 μg
IDH2 (R140Q), FLAG-tag	71100-2	100 μg
IDH1, FLAG-tag	71075-1	50 μg
IDH1, FLAG-tag	71075-2	100 μg
IDH1 (R132H), FLAG-tag	71099-1	50 μg
IDH1 (R132H), FLAG-tag	71099-2	100 μg
IDH2(R140Q) Assay Kit	79309	96 rxns.

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: support@bpsbioscience.com