

ab284508 – DPP4 Inhibitor Screening Kit (Fluorometric)

For the screening of potential DPP4 inhibitors.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab284508>

Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Materials Supplied

Item	Quantity	Storage Condition
DPP4 Assay Buffer	25 mL	-20°C
DPP4 Substrate	200 µL	-20°C
DPP4 Enzyme	100 µL	-20°C
DPP4 Inhibitor (Sitagliptin)	50 µL	-20°C

Reagent Preparation

- Before using the kit, spin the tubes prior to opening.

DPP4 Assay Buffer: Warm to room temperature (RT) before use. Store at 4°C or -20°C.

DPP4 Inhibitor: Store at -20°C. Use within 2 months.

DPP4 Enzyme & DPP4 Substrate: Upon thawing aliquot and store at -20°C, use within 2 months.

Assay Protocol

Enzyme Solution preparation:

- For each well, prepare a 50 µL DPP4 Enzyme Solution:

	Enzyme Solution
DPP4 Assay Buffer	49 µL
DPP4 Enzyme	1 µL

Screen test inhibitors, inhibitor control and blank control Preparations:

- Dissolve test inhibitors into proper solvent.
- Dilute to 4X the desired test concentration with DPP4 Assay Buffer.
- For DPP4 Inhibitor Control, dilute 1:9 with DPP4 Assay Buffer

Δ Note: to compare test inhibitors to Inhibitor Control at its IC₅₀, dilute Inhibitor Control 1:99, then use 25 µL).

- Add 25 µL of test inhibitors, DPP4 Inhibitor or DPP4 Assay Buffer into DPP4 enzyme wells as sample screen (S), Inhibitor Control (IC), or Enzyme Control (EC).
- Mix well, and incubate for 10 minutes at 37°C.

Substrate Mix:

- For each well, prepare a total 25 µL Substrate Solution:

	Substrate Solution
DPP4 Assay Buffer	23 µL
DPP4 Substrate	2 µL

- Mix well and incubate at 37°C

Measurement

Measure the fluorescence (Ex/Em = 360/460 nm) in kinetic mode for 15 - 30 minutes at 37°C. Protect from light. Choose two time points (T₁ & T₂) in the linear range of the plot and obtain the corresponding values for the fluorescence (RFU₁ and RFU₂).

Calculation

- Calculate the slope for all samples, including Enzyme Control (EC), by dividing the net ΔRFU (= RFU₂ - RFU₁) values by the time ΔT (= T₂-T₁).
- Calculate % Relative Inhibition as follows:

$$\% \text{ Relative inhibition} = \frac{(\text{slope of EC} - \text{slope of S})}{\text{slope of EC}} \times 100$$

Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.
For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)