

ab283405 – AMCCase Inhibitor Screening Kit (Fluorometric)

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

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab283405>

Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light. Upon opening, use kit within 6 months.

Materials Supplied

Item	Quantity	Storage Condition
AMCCase Assay Buffer	25 mL	-20°C
Acidic Mammalian Chitinase (lyophilized)	1 vial	-20°C
AMCCase Substrate (in DMSO)	25 µl	-20°C
Bisdionin F (5 mM)	20 µl	-20°C

Materials Required, Not Supplied

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

- Multi-well spectrophotometer (fluorescence plate reader)
- 96-well white opaque plate

Reagent Preparation

- Briefly centrifuge small vials prior to opening.

AMCCase Assay Buffer: Warm to 37 °C before use. Store at either 4°C or -20°C.

AMCCase Substrate: Warm to room temperature (RT) before use.

Acidic Mammalian Chitinase (lyophilized): Reconstitute Acidic Mammalian Chitinase in 550 µl AMCCase Assay Buffer and mix thoroughly. Aliquot and store at -20°C. Avoid repeated freeze/thaw. Keep on ice while in use. Use within two months.

Bisdionin F (5 mM): Warm to room temperature (RT) before use.

Assay Protocol

- **Test Compound preparation:**
 1. Dissolve test sample to 100X in a proper solvent.
 2. Further dilute to 10X in AMCCase Assay Buffer.
 3. Prepare a 20-fold dilution of Bisdionin F i.e. Add 2 µl of the 5 mM Bisdionin F stock solution to 38 µl AMCCase Assay Buffer and mix thoroughly.
 4. Add 10 µl diluted test sample or Diluted Bisdionin F or AMCCase Assay Buffer into wells assigned as test sample (sample, S), Inhibitor Control (IC), or AMCCase Enzyme Control (EC) wells, respectively.
 5. Additional wells with serial dilutions of the test sample may be prepared at this time if desired, containing 10 µl in each candidate well.

Δ Note: Various solvents, in which certain inhibitors are dissolved in, can reduce the AMCCase enzyme activity. Prepare parallel well(s) as Solvent Control (SC) to test the effect of the solvent on AMCCase activity. In case SC is significantly different from EC use its values to determine effect of tested compound.

- **AMCCase Enzyme Solution Preparation:**

1. Prepare a 16-fold dilution of Acidic Mammalian Chitinase (i.e. Dilute of 20 µl of Acidic Mammalian Chitinase with 300 µl of AMCCase Assay Buffer), mix thoroughly and keep on ice.
2. Add 80 µl of Diluted AMCCase Enzyme Solution to each well containing test sample, Inhibitor Control, Solvent Control and AMCCase Enzyme Control.
3. Mix well and incubate at 37 °C for 10-15 minutes. Protect from light.

- **AMCCase Substrate Solution Preparation:**

1. Prepare a 45-fold dilution of AMCCase Substrate Stock Solution (i.e. Dilute 1 µl of AMCCase Substrate with 44 µl of AMCCase Assay Buffer), vortex briefly and keep in ice.
2. After incubation time add 10 µl of Diluted AMCCase Substrate Solution to each well containing test sample, Inhibitor Control, Solvent Control and AMCCase Enzyme Control.

Measurement:

Measure fluorescence (Ex/Em = 320/445nm) in kinetic mode at 37 °C for 30 minutes. Choose two points (t_1 and t_2) in the linear range of the plot and obtain the corresponding fluorescence values (RFU₁ and RFU₂).

Calculation:

Calculate the slope for all samples, including Enzyme Control (EC), by dividing the net Δ RFU (RFU₂-RFU₁) values by the time Δt (t_2-t_1). Calculate % Relative Inhibition and % Relative Activity as follows:

$$\% \text{ Relative Inhibition} = \frac{\text{slope of (enzyme control [EC])} - \text{slope of (test sample [S])}}{\text{slope of (enzyme control [EC])}} \times 100$$

$$\% \text{ Relative activity} = \frac{\text{slope of (test sample [S])}}{\text{slope of (enzyme control [EC])}} \times 100$$

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)