

Instruction Manual

Catalog Number	PK-CA577-K160												
Description	Caspases have been shown to play a crucial role in apoptosis induced by various deleterious and physiologic stimuli. Inhibition of caspases can delay apoptosis, implicating a potential role in drug screening efforts. The PromoKine Caspase-10 Inhibitor Drug Screening Kit provides an effective means for screening caspase inhibitors using a fluorometric approach. The assay utilizes a synthetic peptide substrate AEVD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active Caspase-10 cleaves the synthetic substrate to release free AFC which can then be quantified by fluorometry. Compounds to be screened can directly be added to the reaction and the level of inhibition of Caspase-10 activity can be determined by comparison of the fluorescence intensity in samples with and without the testing inhibitors. The assay is simple, straightforward, and can be performed directly in microtiter plates. Each kit contains 100 units of active Caspase-10, sufficient for screening 100 caspase inhibitor samples. Assay conditions have been optimized to obtain the maximal activity.												
Quantity	100 assays												
Kit Components	<table border="1"> <thead> <tr> <th>Components</th> <th>Quantity</th> </tr> </thead> <tbody> <tr> <td>Caspase Substrate AEVD-AFC (1 mM)</td> <td>0.5 ml</td> </tr> <tr> <td>2x Reaction Buffer</td> <td>10 ml</td> </tr> <tr> <td>Active Caspase-10 (lyophilized)</td> <td>100 units</td> </tr> <tr> <td>DTT (1 M)</td> <td>100 μl</td> </tr> <tr> <td>Caspase Inhibitor, Z-VAD-FMK (2 mM)</td> <td>10 μl</td> </tr> </tbody> </table>	Components	Quantity	Caspase Substrate AEVD-AFC (1 mM)	0.5 ml	2x Reaction Buffer	10 ml	Active Caspase-10 (lyophilized)	100 units	DTT (1 M)	100 μ l	Caspase Inhibitor, Z-VAD-FMK (2 mM)	10 μ l
Components	Quantity												
Caspase Substrate AEVD-AFC (1 mM)	0.5 ml												
2x Reaction Buffer	10 ml												
Active Caspase-10 (lyophilized)	100 units												
DTT (1 M)	100 μ l												
Caspase Inhibitor, Z-VAD-FMK (2 mM)	10 μ l												
Applications / Assay Protocol	<p>A. General Considerations & Reagent Preparations</p> <p>After thawing, store the 2X Reaction Buffer at 4°C. Aliquot enough 2X Reaction Buffer for the number of assays to be performed. Add DTT to the 2X Reaction Buffer immediately before use (10 mM final concentration: add 10 μl of 1.0 M DTT stock per 1 ml of 2X Reaction Buffer).</p> <p>Protect AEVD-AFC from light.</p> <p>Reconstitute the Active Caspase-10 in 550 μl 2X Reaction Buffer. Aliquote and immediately store at -70°C.</p> <p>B. Assay Procedure</p> <ol style="list-style-type: none"> 1. Prepare testing sample in dH₂O to a final volume of 50 μl/well. Add 5 μl of Active Caspase-10. Mix well. Prepare a background control by omitting the Active Caspase-10 from the reaction mixture. Prepare a positive inhibition control by adding 1 μl of the Caspase-10 Inhibitor (provided with the kit) instead of your testing inhibitor. 2. Prepare a Master Mix for each assay containing the follows: 45 μl 2X Reaction Buffer (containing 10 mM DTT) 5μl 1 mM AEVD-AFC substrate (50 μM final concentration). 3. Mix well and add 50 μl of the Master Mix to each well to start the reaction. 4. Incubate at 37°C for 0.5-1 hour. 5. Read samples in a fluorescence plate reader equipped with a 400 nm excitation filter and 505 nm emission filter. Comparison of the fluorescence intensity of the testing samples with samples containing no inhibitors to determine the inhibition efficiency of the testing inhibitors. 												
Storage & Stability	Store kit at -20°C (Store Cell Lysis Buffer, 2X Reaction Buffer, and Dilution Buffer at 4°C after opening). All reagents are stable for at least 6 months.												

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.