

## QuantiFluo™ Urokinase Inhibitor Screening Kit (DUKI-100)

### Quantitative Fluorimetric Inhibitor Screening Assay for Urokinase

#### DESCRIPTION

**UROKINASE PLASMINOGEN ACTIVATOR** (*urokinase*, *uPA*) is a key serine protease involved in the degradation of the extracellular matrix that catalyzes the conversion of plasminogen to active plasmin. It acts as a thrombolytic agent to break up blood clots and when over-expressed, has been reported to influence the growth of certain malignant tumors (breast, prostate, etc.). BioAssay Systems' DUKI-100 Kit provides a convenient fluorimetric means to screen for potential urokinase inhibitors. In this assay, the fluorimetric substrate reacts with urokinase and an inhibitor will decrease the fluorescence at  $\lambda_{\text{ex/em}} = 380/450$  nm.

#### KEY FEATURES

**Safe.** Non-radioactive assay.

**Fast.** Assay is completed within a 30 minute reaction time.

**Convenient and high-throughput.** Homogeneous "mix-incubate-measure" type assay. Can be readily automated to assay thousands of samples per day.

#### APPLICATIONS

For evaluation of drugs and screening potential inhibitors of urokinase.

#### KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

**Assay Buffer:** 10 mL                      **10 mM Inhibitor:** 40  $\mu$ L

**Substrate:** 600  $\mu$ L

**Storage conditions.** The kit is shipped at room temperature. Store all components at  $-20^{\circ}\text{C}$  upon receipt. Shelf life: 12 months after receipt.

**Precautions:** Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to the Material Safety Data Sheet for detailed information.

#### PROCEDURES

This assay is based on a kinetic reaction. To ensure identical incubation time, addition of the Working Reagent (*WR*) to samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended. The assay can be run at room temperature. *Note: Enzyme is not included in the kit. An enzyme of interest is required for this inhibitor assay.*

**Reagent Preparation:** Prior to the assay, equilibrate all components to room temperature and briefly centrifuge tubes before opening. The *WR* should be prepared fresh for each assay run.

**Enzyme Preparation:** Enzyme should be prepared in an enzyme buffer, e.g. 50 mM Tris-HCl, pH 7.4, 100 mM NaCl, 0.08% BSA. The following protocol is optimized for Native Human Urokinase Plasminogen Activator from Cell Sciences (Cat # CRU000A). If using a different enzyme, we recommend that you experimentally determine the optimal amount of enzyme to use per well.

**Test Compound Preparation:** Dissolve test compounds in a solvent of choice (e.g. DMSO). It is prudent to first test the tolerance of the solvent by the enzyme of choice. In the example below, human urokinase (Cell Sciences, Cat# CRU000A) was found to tolerate up to 0.3% DMSO. If a reference  $\text{IC}_{50}$  curve is desired, amiloride hydrochloride inhibitor can be prepared as a test compound according to the concentrations found in the example  $\text{IC}_{50}$  curve.

#### Inhibitor Screening in 96-Well Plate

- Transfer 5  $\mu$ L of Enzyme into separate wells of a black, flat-bottom 96-well plate. Reserve at least one well for No Inhibitor Control ("Control") and one for the No Enzyme Blank ("Blank"): Add 5  $\mu$ L of Enzyme and 5  $\mu$ L of Assay Buffer to the Control and Blank wells respectively.
- To the Control and Blank wells, add 25  $\mu$ L of the solvent in Assay Buffer that the test compounds are dissolved in. For example, if the test compounds are dissolved in Assay Buffer containing 0.1% DMSO, add 25  $\mu$ L of this solution to these wells.

- To the remainder of the wells containing Enzyme, add 25  $\mu$ L of the test compounds. Tap plate to mix and incubate for 10 min at RT to allow the inhibitor to block Enzyme activity.
- Prepare enough *WR* for all wells by mixing 5  $\mu$ L of Substrate and 95  $\mu$ L of Assay Buffer for each well. Transfer 90  $\mu$ L of *WR* to all wells. Briefly tap plate to mix. Incubate for 30 min at room temperature.
- Read the fluorescence intensity at  $\lambda_{\text{ex/em}} = 380/450$  nm.

#### CALCULATION

The percent of Urokinase activity in the presence of test compounds is calculated as follows:

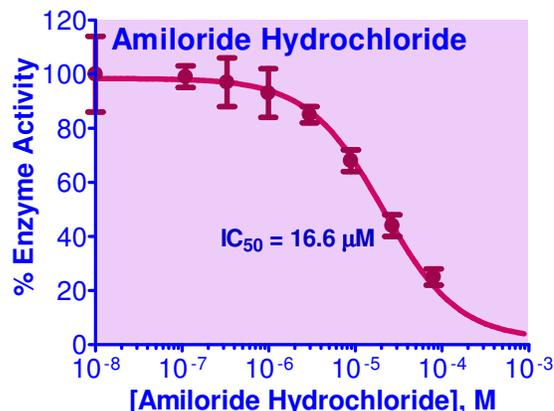
$$\text{Enzyme Activity \%} = \frac{(F_{\text{Compound}} - F_{\text{Blank}})}{(F_{\text{Control}} - F_{\text{Blank}})} \times 100\%$$

Where  $F_{\text{Compound}}$ ,  $F_{\text{Control}}$  and  $F_{\text{Blank}}$  are the fluorescence values of the test compound, Control, and Blank wells at 30 min.

**Inhibitor Screening in 384-Well Plate:** The procedure is essentially the same as for the 96-well plate assay, except that 5  $\mu$ L Enzyme and 15  $\mu$ L of Test Compounds are incubated for 10 min, then 60  $\mu$ L *WR* is added.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices and accessories (e.g. multi-channel pipettor), black, flat-bottom 96-well plates (e.g. Corning Costar), centrifuge tubes and a plate reader.



Inhibition of human urokinase activity by amiloride. Assays are performed in duplicate in 96-well plate according to the standard protocol. Enzyme was incubated with amiloride in the presence of 0.3% DMSO.

#### LITERATURE

- Law, B. et al. (2004). Design, Synthesis, and Characterization of Urokinase Plasminogen-Activator-Sensitive Near-Infrared Reporter. *Chemistry & Biology*. 11, 99–106.
- Mahmood N, Mihalciou C and Rabbani SA (2018) Multifaceted Role of the Urokinase-Type Plasminogen Activator (uPA) and Its Receptor (uPAR): Diagnostic, Prognostic, and Therapeutic Applications. *Front. Oncol.* 8, 24.
- Rockway TW, Nienaber V, Giranda VL. (2002). Inhibitors of the protease domain of urokinase-type plasminogen activator. *Curr Pharm Des.* 8(28), 2541-58.

