

## Amplite™ Fluorimetric Neuraminidase Assay Kit

### \*Blue Fluorescence\*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 12602 (200 assays)	Keep at -20 °C Avoid moisture and light	Fluorescence microplate readers

### Introduction

Neuraminidases, also called sialidases, are glycoside hydrolase enzymes that catalyze the hydrolysis of terminal sialic acid residues and neuraminic acids. The most commonly known neuraminidase is the viral neuraminidase. The cleavage of linkage between sialic acid and adjacent sugar residue permits the transport of the virus through mucin and destroys the haemagglutinin receptor on the host cell, thus allowing elution of progeny virus particles from infected cells. Neuraminidase promotes influenza virus release from infected cells and facilitates virus spread within the respiratory tract. Thus, it is an important target for influenza drug development. The detection of neuraminidase and screening its inhibitors is one of the essential tasks for investigating biological processes and prevention of influenza infection. There are a few assay kits available for detecting neuraminidase, but all the commercial available kits are tedious to use.

Our Amplite™ Fluorimetric Neuraminidase Assay Kit provides a sensitive and robust fluorimetric assay to detect neuraminidase that exists either in cells or biological samples. The non-fluorescent neuraminidase substrate becomes strongly fluorescent upon neuraminidase cleavage. The kit can detect as little as 0.3 mU/mL neuraminidase in a 100 µL assay volume. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. The signal can be easily read by a fluorescence microplate reader at Ex/Em = ~320/~450 nm.

#### Kit Key Features

<b>Broad Application:</b>	Can be used for quantifying neuraminidase in a variety of biological fluids and cells.
<b>Sensitive:</b>	Detect as low as 0.3 mU/mL neuraminidase.
<b>Continuous:</b>	Easily adapted to automation without a separation step.

### Kit Components

Components	Amount
Component A: FluLite™ Blue	1 vial
Component B: Assay Buffer	1 bottle (20 mL)
Component C: Neuraminidase Standard	1 vial (0.1 U)

### Assay Protocol for One 96-well Plate

#### Brief Summary

**Prepare FluLite™ Blue reaction mixture (50 µL) → Add neuraminidase standards or test samples (50 µL)  
→ Incubate at 37 °C or room temperature for 1-2 hours → Monitor fluorescence  
increase at Ex/Em = ~320/~450 nm**

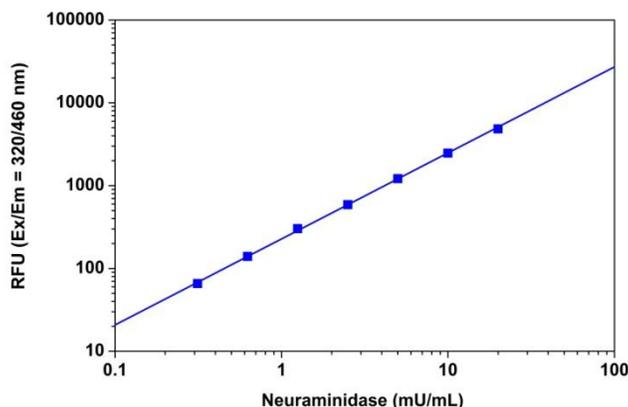
*Note: Thaw all the kit components at room temperature before starting the experiment.*



## Data Analysis

The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the neuraminidase reactions. A neuraminidase standard curve is shown in Figure 1.

*Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*



**Figure 1.** Neuraminidase dose response was measured in a 96-well black plate with Amplite™ Fluorimetric Neuraminidase Assay Kit using a Gemini fluorescence microplate reader (Molecular Devices). As low as 0.3 mU/mL of neuraminidase can be detected with 1 hour incubation time in 37°C, 5% CO<sub>2</sub> incubator.

## References

1. FA Quioco, Carbohydrate-binding proteins: tertiary structures and protein-sugar interactions, Annual Review of Biochemistry, 1986, Vol. 55: 287-315, 1986
2. W. G. Laver, P. M. Colman, R. G. Webster, V. S. Hinshaw and G. M. Air, Influenza virus neuraminidase with hemagglutinin activity, Virology, Vol.137, ( 2), 314-323
3. JN Varghese, WG Laver, PM Colman, Structure of the influenza virus glycoprotein antigen neuraminidase at 2.9 Å resolution, Nature, 303, 35-40, 1983
4. PM Colman, JN Varghese, WG Laver, Structure of the catalytic and antigenic sites in influenza virus neuraminidase, Nature, 303, 41-44, 1983
5. PM Colman, WG Laver, JN Varghese, AT Baker, Three-dimensional structure of a complex of antibody with influenza virus neuraminidase, Nature, 326, 41-44, 1987

**Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.**