

NA-Fluor™ Influenza Neuraminidase Assay Kit: Standardization of the classic MUNANA-based neuraminidase activity assay for monitoring antiviral resistance in influenza



Albana Mihali, Brian D'Eon, and Melissa A. Gee*, Life Technologies, Inc., 35 Wiggins Avenue, Bedford, MA 01730 U.S.A.

ABSTRACT

Here, we describe a MUNANA-based assay kit, the NA-Fluor™ Influenza Neuraminidase Assay Kit, which provides standardization and reproducibility of IC₅₀ methodology. The assay utilizes MUNANA substrate concentration and other conditions comparable to protocols that have been used historically to generate influenza antiviral resistance data. We have benchmarked this assay against the four different MUNANA-based protocols provided on the Neuraminidase Inhibitor Susceptibility Network (NISN) website, http://www.nisn.org/au_about_us.html. The NA-Fluor™ assay gives similar IC₅₀ values for oseltamivir carboxylate and zanamivir when compared to NISN protocols for several tested oseltamivir sensitive and resistant viral strains and can be used to detect mutant virus in mixed viral populations. The assay can be performed in either the traditional endpoint format or using a real-time, kinetic read and is compatible with typical methods to reduce viral contamination such as the addition of detergents or ethanol. The assay is easy to use, reproducible with a Z' of 0.8, and the fluorescence signal is stable for hours making this assay compatible with processing many different viral isolates in a surveillance screen or for high throughput screening for drug lead identification.

INTRODUCTION

Although neuraminidase inhibitors (NIs) remain an effective treatment for influenza virus strains currently considered world health threats, such as Influenza virus A which includes pandemic H1N1 and Influenza virus B, viral resistance to these drugs threatens their usefulness for future pandemics. Key antivirals oseltamivir carboxylate (Tamiflu®) and zanamivir (Relenza®) target the viral enzyme neuraminidase which is responsible for release of newly generated viral particles from an infected cell's surface, a process necessary for transmission of the virus to other cells within the organism and for transmission between organisms. Oseltamivir resistance in seasonal H1N1 increased from 12.3% during the 2007-2008 flu season to 98.5% in 2008-2009 season (1), heightening concern that drug resistance will likewise become prominent in pandemic viral strains and highlighting the need for antiviral drug resistance surveillance. The H275Y mutation in H1N1 neuraminidase is the most common mutation conferring resistance, however, due to the high mutation rates of viruses, new mutations can be expected that will also render viral neuraminidase less sensitive to antiviral drugs. PCR methods can be used to detect previously identified mutations, however, functional neuraminidase enzyme activity inhibition testing is necessary for detecting drug resistance that results from novel mutations. The two neuraminidase enzyme inhibition assays using either the fluorescent MUNANA or chemiluminescent NA-Star® substrate are robust tools for NI susceptibility testing. The MUNANA-based assay is broadly used by many groups including several world health organizations for NI susceptibility testing, yet no standardized protocol or dedicated kit is in place for this assay, making comparison of data generated between different laboratories difficult.

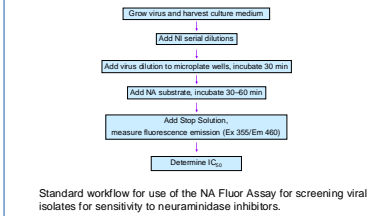
MATERIALS AND METHODS

All assays, unless otherwise noted, were conducted using the standard NA-Fluor™ Influenza Neuraminidase Assay Kit protocol. Basically, NIs were prepared in 10-fold serial dilutions at 4x final concentration in assay buffer (16.65 mM MES, 2mM CaCl₂, pH 6.5) and 25 µl were added to wells of a black flat-bottom 96-well microplate. Viral samples were diluted (1:3 to 1:100 depending on virus strain and titer) in assay buffer and 25 µl were added to the NI serial dilutions for preincubation (30 min at 37°C). NA-Fluor™ Substrate was added at 50 µl for a final assay concentration of 100 µM MUNANA and incubated (37°C for 60 min). No-virus controls were included on each assay plate. Assay was terminated by addition of 100 µl of NA-Fluor™ Stop Solution and plates were read on a SpectraMax® M5 or M2 (Molecular Devices) plate reader using Ex 355nm/Em 460nm settings. For data analysis, no-virus control assay well RFLUs were subtracted from viral data RFLUs and data was processed using either Microsoft Excel or GraphPad® Prism software. For real-time detection, the assay was performed as above but without the stop solution addition and assay reads were taken at 5 min intervals.

Viral samples were titrated according to the NA-Fluor™ Assay protocol. Influenza A/WS/33 (H1N1) (VR-1520™) and influenza B/Lee/40 (VR-1535™) were purchased from ATCC. Other viral strains were gifts (See Acknowledgements.)

RESULTS

Figure 1. Neuraminidase Inhibition Assay Workflow



Comparison to Traditional MUNANA Assays

Figure 2. Comparison of NA-Fluor Assay Sensitivity

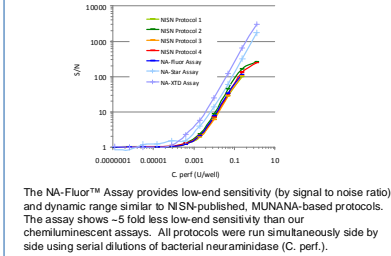
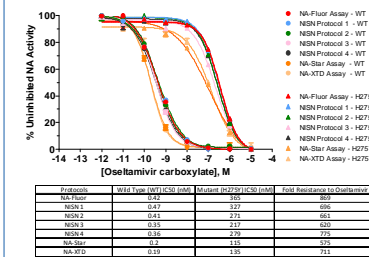
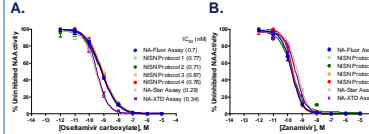


Figure 3. Comparison of NA-Fluor™ Assay to Other Assays for IC₅₀ Determination



The NA-Fluor™ Assay provides IC₅₀ values and sensitivity for detecting sensitive virus similar to NISN-published, MUNANA-based protocols. All protocols were run in parallel to determine the IC₅₀ values of influenza A/Texas/36/91 (H1N1) sensitive and oseltamivir-resistant mutant (H275Y) strains for oseltamivir carboxylate.

Figure 4. Additional Protocol Comparison



Compatibility for High Throughput Processing

Figure 5. NA-Fluor™ Assay Signal Stability with Time

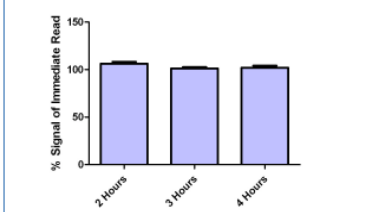


Table 1. Reproducibility of NA-Fluor™ Assay

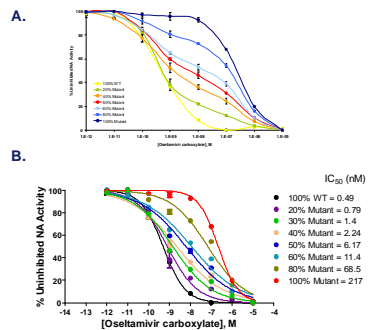
Time (hours)	0	2	3	4
NA-Fluor Assay % Read-out (96 reps)	100	106.3	101.2	102.0
St. Dev. (16 no virus reps, 40 no drug reps, 40 10nM drug reps)	0	6.0	1.4	2.2
Z' = 1 - (3σ ₁ + 3σ ₂) / (μ ₁ - μ ₂)	0.78	0.80	0.80	0.80

σ = St. Dev.
μ = mean
σ₁ = no inhibitor
σ₂ = 100% inhibitor

The NA-Fluor™ assay is highly reproducible giving a Z' of 0.78 or above indicating the assay can be used confidently to identify neuraminidase inhibitors in high throughput screening mode.

Detection of Mutants in Mixed Viral Populations

Figure 6. Detection of NI-Resistant H1N1 Virus in Mixed Populations



Assay Flexibility

Figure 7. Assay Time Flexibility

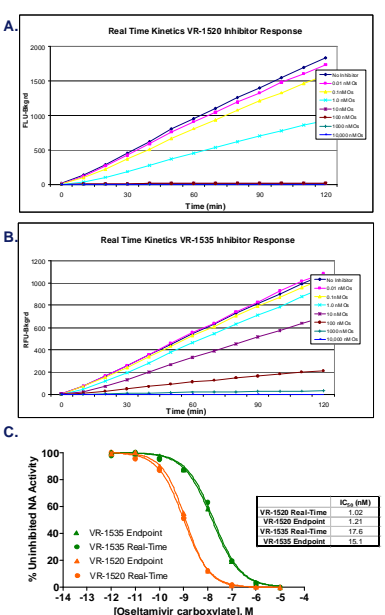


Figure 8. NA-Fluor™ Assay Flexibility in Time and Virus Dilution

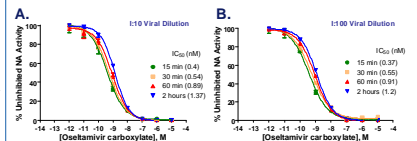
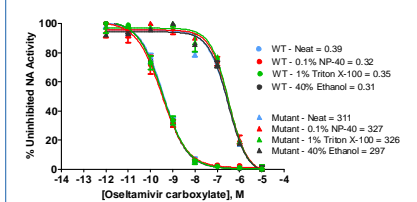


Figure 9. Compatibility with Viral Inactivation Methods.



CONCLUSIONS

We have developed a standardized MUNANA-based neuraminidase assay kit, the NA-Fluor™ Influenza Neuraminidase Assay Kit, that has been optimized for NI susceptibility screening. The assay is:

- A standardized MUNANA-based fluorescent assay.
- Economically priced
- Provides data that can be compared to data generated using traditional MUNANA-based protocols
- Has stable signal read-out allowing for high throughput processing
- Easy to use
- Highly reproducible
- Provides flexibility in assay format
- Compatible with viral inactivation methods
- Environmentally friendly

REFERENCES

1. Dharan, NU et al. 2009. J American Medical Association. 301:1034-1041.
2. NISN IC50 Methodology (<http://www.nisn.org/>); Marathe, BM, and EA Govorkova (2010). Determination of influenza virus susceptibility to neuraminidase inhibitors using fluorescent substrate. St. Jude Children's Research Hospital.
3. Hurt, A. (2009) Fluorometric neuraminidase inhibition assay. Standard Operating Procedure. WHO-025. WHO Collaborating Centre for Reference and Research on Influenza, Australia.
4. McKimm-Breschkin, JA. Enzyme inhibition assay for sensitivity to NI inhibitors. Fluorescence based assay (MUNANA). Lackerby, A. (2008). Determination of influenza virus susceptibility to neuraminidase inhibitors using a fluorescent substrate.
5. Zhang JH, et al. 1999. J Biomedical Screening. 4:67-73.
6. Jonges, M., et al. 2010. J Clinical Microbiology. 48:928-940.

ACKNOWLEDGEMENTS

Influenza A/Texas/36/91 (H1N1) wild-type and H274Y strains were kindly provided by Dr. Larisa Gubareva (Influenza Branch, CDC, USA). Oseltamivir carboxylate was kindly provided by Hoffman-La Roche Inc.. Zanamivir was kindly provided by GlaxoSmithKline.

TRADEMARKS/LICENSING

Applied Biosystems, AB Design, Invitrogen, Life Technologies, NA-Fluor, NA-Star and NA-XTD are trademarks of Life Technologies. All other trademarks are the property of their respective owners.

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.