

ab185436

**Neuraminidase Activity
Assay Kit (Fluorometric)**

Instructions for Use

For the rapid, sensitive and accurate measurement of Neuraminidase activity in various samples.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

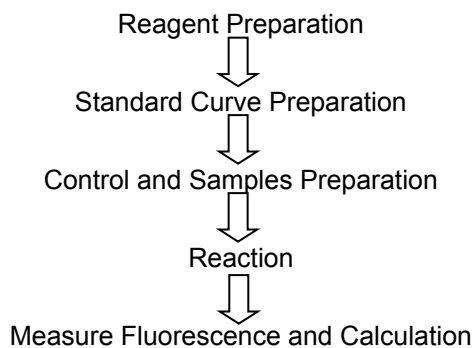
1.	Overview	3
2.	Protocol Summary	4
3.	Kits Components	5
4.	Storage and Stability	6
5.	Materials Required, Not Supplied	6
6.	Reagents Preparation	7
7.	Assay Protocol	8
8.	Data Analysis	11
9.	Troubleshooting	15

1. Overview

Neuraminidase (NA) is a very common enzyme that hydrolyzes terminal sialic acid residues on polysaccharide chains; most often a galactose residue. NA activity plays a key role in the invasion of target cells and the replication of influenza virus. NA activity also assists in the elution of progeny viruses from infected cells, and prevents self-aggregation of virus. Thus, NA is an important target for drug development.

Abcam's Neuraminidase Activity Assay Kit (Fluorometric) (ab185436) provides a simple and sensitive method for measuring NA activity using fluorescence (Ex/Em = 530/590 nm). The assay utilizes NA Probe to detect the neuraminidase activity. This high-throughput adaptable assay kit can detect NA activity as low as 2.0 mU/ml in a variety of samples.

2. Protocol Summary



3. Kits Components

Item	Quantity
NA Assay Buffer	30 mL
NA Probe	200 µL
NA Enzyme Mix I	1 vial
NA Enzyme Mix II	1 vial
Galactose Standard (100 nmol/µL)	100 µL
NA Positive Control	100 µL
NA Substrate	1 vial

4. Storage and Stability

Upon arrival, store the kit at -20°C and protect from light. Please read the entire protocol before performing the assay. Avoid repeated freeze/thaw cycles.

Briefly centrifuge all small vials prior to opening.

5. Materials Required, Not Supplied

- Distilled water (dH₂O)
- 96-well black plate with flat bottom
- Multi-well spectrophotometer (ELISA reader)

6. Reagents Preparation

1. NA Assay Buffer and NA Probe:

Warm to room temperature before use. Aliquot and store at -20°C. Stable for two months.

2. NA Substrate:

Reconstitute with 110 µL dH₂O. Aliquot and store at -20°C. Stable for two months.

3. NA Positive Control:

Aliquot and store at -20°C. Use within two months

4. NA Enzyme Mix I and Enzyme Mix II:

Reconstitute with 220 µL NA Assay Buffer. Aliquot and store at -20°C. Stable for two months.

7. Assay Protocol

1. Sample Preparation

- a) Homogenize 10 mg of sample (wet weight or cell pellet) in 100 μ L NA Assay Buffer.
- b) Centrifuge at 10,000 x *g* for 5 minutes at 4°C. Collect the supernatant.
- c) Add 2-10 μ L of supernatant or serum into a 96-well plate and adjust the volume to 50 μ L with NA Assay Buffer.
- d) Add 10 μ L of NA Positive Control into desired wells(s) and adjust the volume to 50 μ L with NA Assay Buffer.

NOTE:

- *For unknown samples, we suggest performing pilot experiment & testing several doses of your samples to ensure the readings are within the Standard Curve range.*
- *For samples having background, prepare parallel well(s) containing the same amount of sample as in the test well (sample background control). Adjust the volume to 50 μ L with NA Assay Buffer.*
- *1% or higher Triton X-100 concentration interferes with the assay.*

2. Standard Curve Preparation:

- a) Dilute Galactose Standard to 1 nmol/ μ L by adding 10 μ L of 100 nmol/ μ L Galactose Standard into 990 μ L of NA Assay Buffer, mix well. Dilute the Standard further to 0.1 nmol/ μ L by adding 20 μ L of 1 nmol/ μ L Standard to 180 μ L of NA Assay Buffer and mix.
- b) Add 0, 1, 2, 3, 4 and 5 μ L of 0.1 nmol/ μ L Galactose Standard into a series of wells in 96-well plate to generate 0, 0.1, 0.2, 0.3, 0.4 and 0.5 nmol/well of Galactose Standard.
- c) Adjust the volume to 50 μ L/well with NA Assay Buffer.

3. Reaction Mix:

Prepare enough Reaction Mix for the number of assays to be performed. For each well, prepare 50 μ L Mix containing:

	Reaction Mix	Background Control Mix
NA Assay Buffer	44.5 μ L	45.5 μ L
NA Enzyme Mix I	2 μ L	2 μ L
NA Enzyme Mix II	2 μ L	2 μ L
NA Substrate	1 μ L	-----
NA Probe	0.5 μ L	0.5 μ L

Add 50 μL of Reaction Mix to each well containing the Standards, Positive Control and samples. Mix well.

* For samples having background, add 50 μL of Background Control Mix to sample background control well(s). Mix well.

4. Measurement

Incubate for 30 minutes at 37°C and measure fluorescence (Ex/Em = 535/590 nm) kinetically.

NOTE:

- *Incubation time depends on the NA Activity in the samples. Choose two time points (T1 and T2) in the linear range (fluorescence values A1 and A2 respectively) to calculate the NA activity of the samples. The Standard Curve can be read in end point mode (i.e. at the end of incubation time).*

8. Data Analysis

Calculations:

- a) Subtract 0 Standard reading from all readings. Plot the Galactose Standard Curve. If sample background control reading is significant, subtract background control reading from sample readings. Calculate the NA activity of the test sample: $\Delta\text{RFU} = A_2 - A_1$. Apply ΔRFU to the Standard Curve to get B nmol of Galactose generated by NA during the reaction time ($\Delta T = T_2 - T_1$).

$$\text{Sample Neuraminidase Activity} = \frac{B}{(\Delta T \times V)} \times D = \text{nmol/min/ml} \\ = \text{mU/ml}$$

Where:

B is the Galactose amount from the Standard Curve (nmol)

ΔT is the reaction time (minutes)

V is the sample volume added into the reaction well (mL)

D is the sample dilution factor

Unit Definition: One unit of Neuraminidase activity is the amount of enzyme that generates 1.0 μmol of Galactose per min. at pH 7.4 at 37°C.

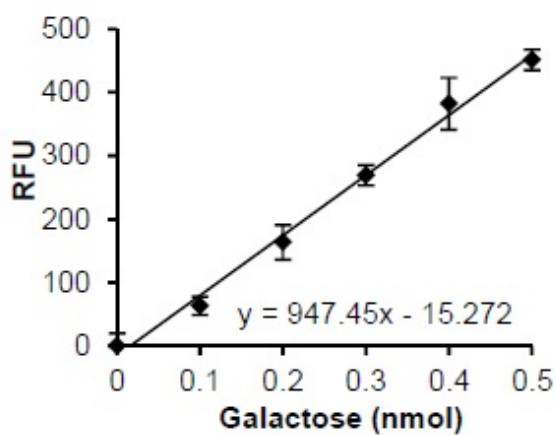


Figure 1. Galactose Standard Curve.

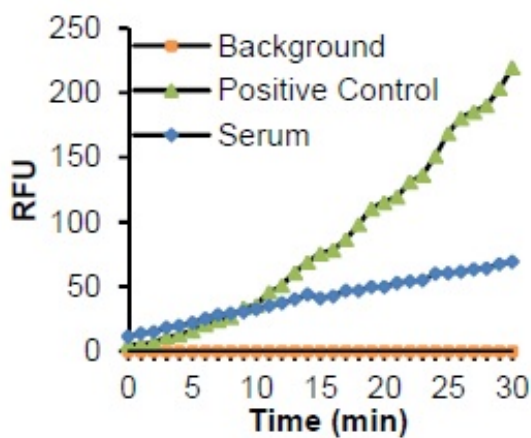


Figure 2. NA activity in normal Human serum (1 μ L) and positive control (1 μ L). Assays were performed following the kit protocol.

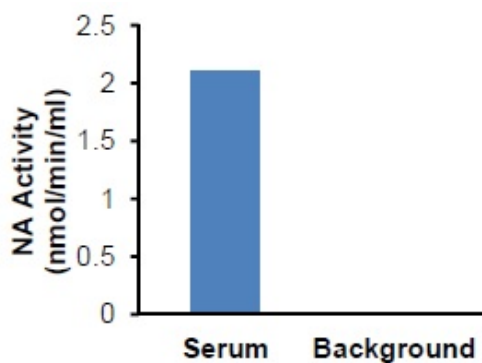


Figure 3. Calculated activity of serum. Assays were performed following the kit protocol.

9. Troubleshooting

Problem	Reason	Solution
Assay not working	Assay buffer at wrong temperature	Assay buffer must not be chilled - needs to be at RT
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended samples types as listed on the datasheet
	Sample readings are outside linear range	Concentrate/ dilute samples to be in linear range

Problem	Reason	Solution
Samples with inconsistent readings	Unsuitable sample type	Refer to datasheet for details about incompatible samples
	Samples prepared in the wrong buffer	Use the assay buffer provided (or refer to datasheet for instructions)
	Samples not deproteinized (if indicated on datasheet)	Use the 10kDa spin column (ab93349)
	Cell/ tissue samples not sufficiently homogenized	Increase sonication time/ number of strokes with the Dounce homogenizer
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Samples are too old or incorrectly stored	Use freshly made samples and store at recommended temperature until use
Lower/ Higher readings in samples and standards	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Out-of-date kit or incorrectly stored reagents	Always check expiry date and store kit components as recommended on the datasheet
	Reagents sitting for extended periods on ice	Try to prepare a fresh reaction mix prior to each use
	Incorrect incubation time/ temperature	Refer to datasheet for recommended incubation time and/ or temperature
	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)

Problem	Reason	Solution
Standard curve is not linear	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Pipetting errors when setting up the standard curve	Try not to pipette too small volumes
	Incorrect pipetting when preparing the reaction mix	Always prepare a master mix
	Air bubbles in wells	Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates
	Concentration of standard stock incorrect	Recheck datasheet for recommended concentrations of standard stocks
	Errors in standard curve calculations	Refer to datasheet and re-check the calculations
	Use of other reagents than those provided with the kit	Use fresh components from the same kit

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 108008523689 (中國聯通)

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp