

Product datasheet

Sialic Acid (NANA) Assay Kit ab83375

★★★★★ [1 Abreviews](#) [9 References](#) [3 Images](#)

Overview

Product name	Sialic Acid (NANA) Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Urine, Cell culture media
Assay type	Quantitative
Sensitivity	1 μ M
Range	0.1 μ M - 10 μ M
Assay time	0h 40m
Product overview	Sialic Acid (NANA) Assay Kit (ab83375) is a simple and convenient kit for measuring free sialic acid (mainly N-acetylneuraminic acid (NANA)) in biological samples.

The sialic acid assay uses an enzyme coupled reaction in which oxidation of free sialic acid creates an intermediate that reacts stoichiometrically with the probe to generate a product that can be detected by absorbance (OD = 570 nm) or fluorescence (Ex/Em=535/587 nm).

The kit measures sialic acid in the linear range of 0.1 to 10 nmol with a detection sensitivity ~1 μ M concentration.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K566 Sialic Acid (NANA) Colorimetric/Fluorometric Assay Kit. K566-100 is the same size as the 100 test size of ab83375.

Sialic acid is a generic term for the N- or O-substituted derivatives of neuraminic acid, a monosaccharide with a nine-carbon backbone. It is also the name for the most common member of this group, N-acetylneuraminic acid (NANA). Sialic acids are found widely distributed in animal tissues and to a lesser extent in other species ranging from plants and fungi to yeasts and bacteria, mostly in glycoproteins and gangliosides. It has been shown recently that sialic acid level may be associated with developmental and pathological stages.

Platform Microplate reader

Properties

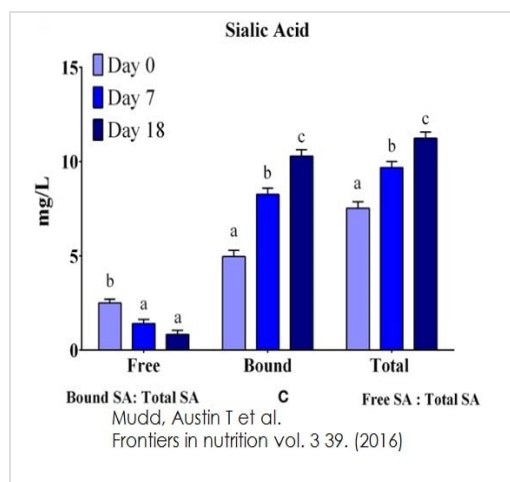
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer IV	1 x 25ml
Development Enzyme Mix I	1 vial
OxiRed Probe	1 x 200µl
Sialic Acid Assay Buffer	1 x 25ml
Sialic Acid Converting Enzyme	1 vial
Sialic Acid Standard	1 vial

Relevance

Sialic acid is a generic term for the N- or O-substituted derivatives of neuraminic acid, a monosaccharide with a nine-carbon backbone. It is also the name for the most common member of this group, N-acetylneuraminic acid (NANA). Sialic acids are found widely distributed in animal tissues and to a lesser extent in other species ranging from plants and fungi to yeasts and bacteria, mostly in glycoproteins and gangliosides. It has been show recently that sialic acid level may associate with developmental and pathological stages.

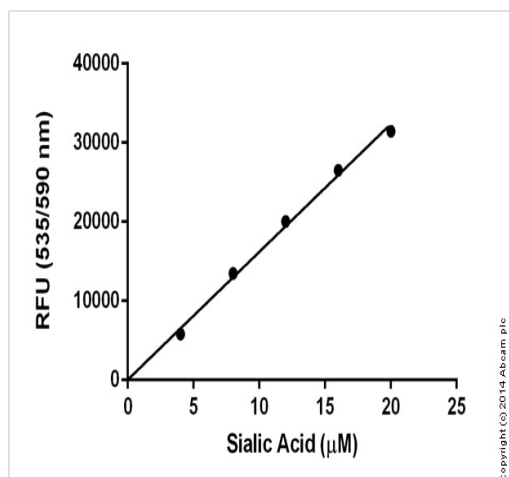
Images



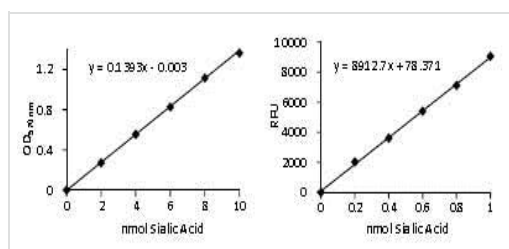
Concentrations of free SA decreased ($P < 0.0001$), whereas bound and total SA increased ($P < 0.0001$) across lactation. Measured using ab83375.

Functional Studies - Sialic Acid (NANA) Assay Kit
(ab83375)

Mudd, Austin T et al., Frontiers in nutrition?vol. 3 39., Fig 3, doi:10.3389/fnut.2016.00039



Functional studies - Sialic Acid (NANA) Assay Kit
(ab83375)



Functional Studies - Sialic Acid (NANA) Assay Kit
(ab83375)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise,

please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors