

## Product datasheet

### 100X NAD+ ab252275

#### Overview

<b>Product name</b>	100X NAD+
<b>Product overview</b>	100X NAD+ is available as a stand-alone component for our Mitochondrial Aldehyde Dehydrogenase (ALDH2) Activity Assay Kit ( <a href="#">ab115348</a> ).  Rehydrate by adding 250 µL deionized water.
<b>Tested applications</b>	<b>Suitable for:</b> EIA

#### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 units
100X NAD+	1 unit

#### Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab252275 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
EIA		Use at an assay dependent concentration.

**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