

Product datasheet

7-AAD Staining Solution ab228563

Overview

Product name	7-AAD Staining Solution
Sample type	Suspension cells
Assay type	Cell-based
Product overview	<p>7-Amino-Actinomycin D (7-AAD) is ready-to-use nucleic acid dye solution (2 ml). 7-AAD can be used in place of propidium iodide (PI) for the exclusion of non-viable cells in flow cytometry analysis. This product can be used in combination with PE (phycoerythrin), and FITC conjugated antibodies in 2-color analysis. The advantage of 7-AAD over PI is the minimal spectral overlap between these emissions. Fluorescence is detected in the far red range of the spectrum (650 nm long-pass filter).</p> <p>Excitation = 540 nm / emission = 645 nm.</p>
Tested applications	Suitable for: Flow Cyt

Properties

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

Components	400 tests
7-AAD Staining Solution	1 x 2ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab228563 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
Flow Cyt		Use 5µl for 10 ⁶ cells. Incubate for 5 minutes at room temperature before analysis.

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