invitrogen by Thermo Fisher Scientific

eBioscience™ 7-AAD Viability Staining Solution

Catalog Number: 00-6993 Also known as: 7AAD, 7-amino-actinomycin D For Research Use Only. Not for use in diagnostic procedures.

Product Information

Contents: eBioscience™ 7-AAD Viability Staining Solution REF Catalog Number: 00-6993

Formulation: PBS, 0.09% sodium azide Temperature Limitation: Store at 2-8°C. Light sensitive material. Batch Code: Refer to vial

- Use Bv: Refer to vial
- Caution, contains Azide

Description

7-AAD (7-amino-actinomycin D) is a ready-to-use solution for the exclusion of nonviable cells in flow cytometric analysis. 7-AAD can be used in place of PI (propidium iodide). This solution can be used in combination with PE (phycoerythrin) and FITC (fluorescein isothiocyanate) conjugated antibodies in 2-color analysis. The advantage of 7-AAD over PI is that there is minimal spectral overlap between these emissions. Fluorescence is detected in the far red range of the spectrum (650 nm long-pass filter).

X

LOT

Applications Reported

7-AAD can be used as a viability probe for methods of nonviable cell exclusion, based on light scatter and uptake of the reagent as detected in FL3. CAUTION: 7-AAD is a potential carcinogen. It is recommended that the user wear protective clothing, gloves, and eye/face protection in order to avoid contact with skin and eyes.

Applications Tested

This lot of 7-AAD has been tested by flow cytometric analysis. Use at 5 μ L (0.25 μ g) per test (per million cells) and incubate for 5 minutes before analysis. A test is defined as the amount (μ g) of antibody that will stain a cell sample in a final volume of 100 μ L. Cell number should be determined empirically but can range from 10⁵ to 10⁸ cells/test.

References

MC O'Brien, et al. (1995) Cytometry 19:243 I Schmid et al. (1992) Cytometry 13:204

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com