

7-AAD

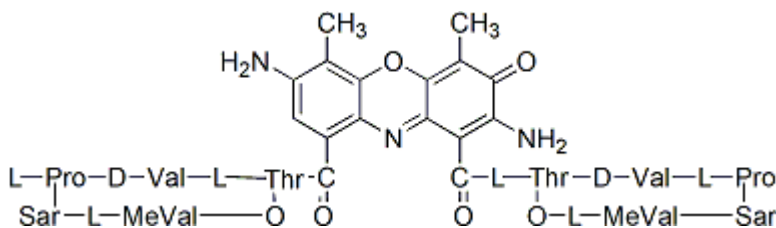
7-AAD (7-aminoactinomycin D) is a membrane-impermeant fluorescent DNA binding dye that is useful for live/dead discrimination and cell cycle profiling by flow cytometry.



Product Description

7-AAD is a fluorescent DNA binding dye that is membrane impermeant and therefore generally excluded from live cells and early apoptotic cells, but stains necrotic and late apoptotic cells with compromised membrane integrity.

- Selective detection of dead cells by flow cytometry in the PE-Cy5/PerCP channel
- Perform cell cycle profiling by flow cytometry in fixed/permeabilized cells
- $\lambda_{Ex}/\lambda_{Em}$ (with DNA) = 546/647 nm (7-AAD can also be excited by the 488 or 532 laser lines)
- Orange red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light, especially in solution
- $C_{62}H_{87}N_{13}O_{16}$
- MW: 1270.45



7-AAD bound to DNA has a large Stokes shift with excitation/emission at 546/647 nm. The dye can be excited with either the 488 nm or 532 nm laser lines, and detected in the PE-Cy5/PerCP flow cytometry channel. 7-AAD is widely used for live/dead cell discrimination. It also can be used in fixed and permeabilized cells for cell cycle profiling by DNA content analysis using flow cytometry. 7-AAD intercalates selectively at GC-rich regions of DNA, making it dye useful for chromosome banding studies. We offer 7-AAD as a solid powder (catalog no. 40037) or as a solution in 1 mg/mL DMSO:Water (1:1) (catalog no. 40084).

7-AAD products

Product	Catalog Number	Unit Size	Format
7-AAD	40037	1 mg	Orange/red solid
7-AAD Solution, 1 mg/mL	40084	1 mL	Red solution in DMSO:Water (1:1)

We also offer the [CF@488A Annexin V and 7-AAD Apoptosis kit](#), containing CF@488A Annexin V for staining apoptotic cells green, and 7-AAD for staining necrotic cells with far-red fluorescence. Also see [NucSpot@ Far-Red](#), an improved alternative to 7-AAD that has red-shifted fluorescence emission, for less bleed-through fluorescence in the PE-Texas Red® channel. For fixable dead cell stains, see our [Live-or-Dye™ Fixable Viability Staining Kits](#).

Cy Dye is a registered trademark of Cytiva; Texas Red is a registered trademark of Thermo Fisher Scientific.

References

1. Environ Mol Mutagen (2013) 54(8), 682-9. [doi: 10.1002/em.21800](https://doi.org/10.1002/em.21800)
2. Mol Cell Biol (2014) 34(24), 4485-4499. [doi: 10.1128/MCB.00671-14](https://doi.org/10.1128/MCB.00671-14)
3. Mutat Res Genet Toxicol Environ Mutagen (2014) 770, 1-5. [doi: 10.1016/j.mrgentox.2014.05.001](https://doi.org/10.1016/j.mrgentox.2014.05.001)
4. Science (2017) 356(6339) eaai9264. [DOI: 10.1126/science.aai9264](https://doi.org/10.1126/science.aai9264)
5. Int J Rad Biol (2019) 95,9, 1226-1235. [DOI: 10.1080/09553002.2019.1625490](https://doi.org/10.1080/09553002.2019.1625490)
6. OncoTargets Therapy (2019) 12, 5227-5239. <https://doi.org/10.2147/OTT.S190460>
7. Antioxidants (2020) 9(9), 873. [doi:10.3390/antiox9090873](https://doi.org/10.3390/antiox9090873)
8. Antioxidants (2020) 9(9), 876. [doi:10.3390/antiox9090876](https://doi.org/10.3390/antiox9090876)

This datasheet was generated on June 20, 2026 at 07:24:33 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/7-aad-7-aminoactinomycin-d/>

Product attributes

Probe cellular localization	Nucleus
Assay type/options	DNA content/cell cycle profiling (flow cytometry), Live/dead discrimination, No-wash staining
Detection method/readout	Flow cytometry
Cell permeability	Membrane impermeant
Apoptosis/viability marker	Dead cell stain
Colors	Far-red
Excitation/Emission	546/647 nm (with DNA)