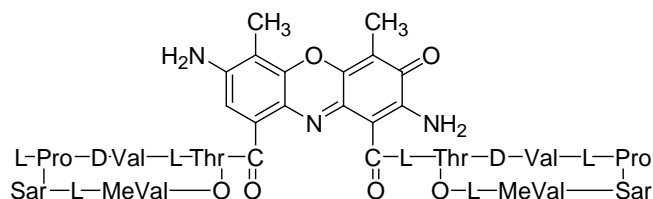


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 7-AAD (7-aminoactinomycin D)**CATALOG #** 40037**MOLECULAR INFORMATION:** C₆₂H₈₇N₁₃O₁₆
MWt: 1270.45**PROPERTIES:**

Color & Form Orange red solid
Purity ≥ 95% by TLC
Solubility Soluble in DMSO or DMF
Absorption/Emission 546/647nm (DNA)
Extinction Coefficient

STORAGE AND HANDLING: Store at -20°C, protect from light.**APPLICATION:** 7-AAD is a fluorescent DNA intercalator, which upon binding to dsDNA can be excited by the argon-ion laser and emits at 647 nm.¹⁻³ 7-AAD binds selectively to GC regions of DNA, which makes the dye useful for chromosome banding studies.⁴

Ref.: 1) *Exp. Parasitol.* **97**, 141 (2001); 2) *Br. J. Haematol.* **104**, 530(1990); 3) *Cytometry* **12**, 221(1991); 4) *Chromosoma* **68**, 287(1978).

TOXICITY: Unknown

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.
