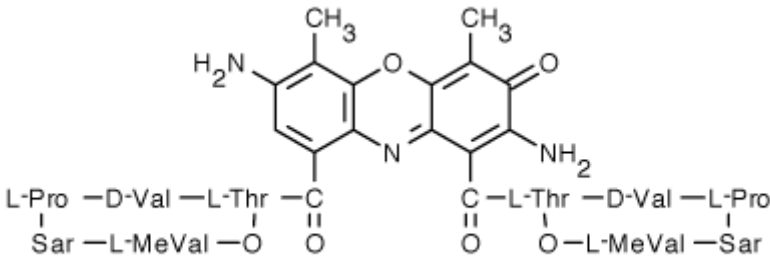


## 7-aminoactinomycin D

## Instruction Manual

Catalog Number	PK-CA707-40037
Description	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. 7-AAD binds selectively to GC regions of DNA, which makes the dye useful for chromosome banding studies.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}(\text{DNA}) = 546/647 \text{ nm}$
Molecular Structure	
Molecular Weight / Molecular Formula	1270.45 Da; $\text{C}_{62}\text{H}_{87}\text{N}_{13}\text{O}_{16}$
Purity	$\geq 95\%$ (as determined by TLC)
Appearance / Formulation / Solubility	Orange red solid; soluble in DMF or DMSO.
Storage & Stability	Store at $-20^{\circ}\text{C}$ and protect from light, especially in solution.
Applications	Cell stain; fluorescent marker for double-stranded DNA (nucleus staining)
References	<ol style="list-style-type: none"> <li>1) Exp. Parasitol. 97, 141 (2001)</li> <li>2) Br. J. Haematol. 104, 530(1990)</li> <li>3) Cytometry 12, 221(1991)</li> <li>4) Chromosoma 68, 287(1978)</li> </ol>
Caution	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.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.