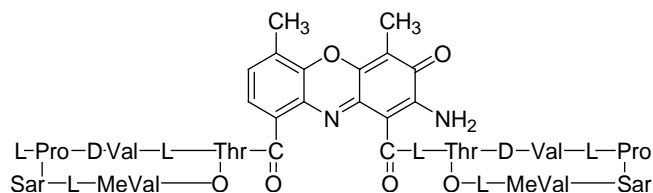


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Actinomycin D**CATALOG #** 40038**MOLECULAR INFORMATION:** C<sub>62</sub>H<sub>86</sub>N<sub>12</sub>O<sub>16</sub>  
MWt: 1255.43**PROPERTIES:**

**Color & Form** Orange red solid  
**Purity** ≥ 98% by HPLC  
**Solubility** Soluble in DMSO at 50mg/mL  
**Absorption/Emission** 240nm & 443nm/none  
**Extinction Coefficient** >35,000(240nm) & >25,000(443nm)

**STORAGE AND HANDLING:** Stored at 4 °C.**APPLICATION:** Actinomycin D is a nonfluorescent GC-selective intercalator.<sup>1</sup> Similar to 7-AAD,<sup>2</sup> actinomycin D has also been used for chromosome banding studies.Ref.: 1) *J. Mol. Biol.* **225**, 445(1992); 2) *Cancer Genet. Cytogenet.* **1**, 187(1980).**TOXICITY:** Unknown

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> 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.
---