

Actinomycin D

Actinomycin D is a nonfluorescent GC-selective intercalator. Similar to 7-AAD (cat. no. 40038), actinomycin D has also been used for chromosome banding studies.



Product attributes

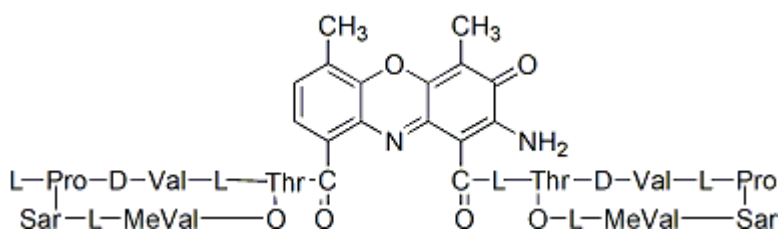
Storage Conditions

Store at 2 to 8 °C

Product Description

Actinomycin D is a nonfluorescent GC-selective intercalator. Similar to 7-AAD (cat. no. 40037), actinomycin D has also been used for chromosome banding studies.

- Orange red solid soluble in DMSO at 50 mg/mL
- Store at 4 °C
- $C_{62}H_{86}N_{12}O_{16}$
- MW: 1255.43



This datasheet was generated on December 5, 2025 at 07:23:21 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/actinomycin-d/>