APTS

APTS is a very useful green fluorescent dye for labeling glycoproteins or sugar molecules in general. The labeling occurs via reductive amination which involves the formation of a Schiff base between the amine of APTS and the aldehyde or ketone of the sugar, followed by the reduction of the Schiff base linkage to a stable carbon-nitrogen bond.

sistum

Product attributes

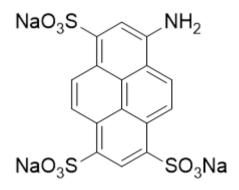
CAS number 196504-57-1 **Excitation/Emission** 424/505 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

APTS is a very useful green fluorescent dye for labeling glycoproteins or sugar molecules in general. The labeling occurs via reductive amination which involves the formation of a Schiff base between the amine of APTS and the aldehyde or ketone of the sugar, followed by the reduction of the Schiff base linkage to a stable carbon-nitrogen bond. The multi-anionic nature of the dye makes it ideal for the studies of carbohydrate molecules by high resolution capillary electrophoresis.

- λ_{Ex}/λ_{Em}: 424/505 nm (before conjugation)
- Yellow solid soluble in water
- Store at 4°C and protect from light.
- C₁₆H₈Na₃O₉S₃
- MW: 523.39
- [196504-57-1]



This datasheet was generated on November 6, 2025 at 03:03:06 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/apts-8-aminopyrene-136-trisulfonic-acid-trisodium-salt/