

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.



Product attributes

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

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.

- $\lambda_{Ex}/\lambda_{Em}$: 424/505 nm (before conjugation)
- Yellow solid soluble in water
- Store at 4 °C and protect from light.
- $C_{16}H_8Na_3O_9S_3$
- MW: 523.39
- [196504-57-1]

