

APTS [8-Aminopyrene-1,3,6-trisulfonic acid, trisodium salt] *CAS 196504-57-1*

Catalog number: 624 Unit size: 10 mg

Component	Storage	Amount
APTS [8-Aminopyrene-1,3,6-trisulfonic acid, trisodium salt] *CAS 196504-57-1*	Freeze (<-15 °C), Minimize light exposure	10 mg

OVERVIEW

This water-soluble reagent has strong green fluorescence with great Stokes shift. It is widely used for detecting carbohydrates and labeling glycoproteins in electrophoresis. It is also used to modify other carbonyl-containing biomolecules such as aldehydes and ketones.

AT A GLANCE

 $\label{eq:monotone} \textbf{Important} \quad \text{It is important to store APTS at <-15 °C and should be stored in cool, dark place.}$

It can be used within 24 months from the receiving date.

KEY PARAMETERS

Instrument: Fluorescence microplate reader

Excitation: 424 nm
Emission: 505 nm
Cutoff: 450 nm
Recommended plate: Solid black

PREPARATION OF WORKING SOLUTION

APTS working solution:

Add water to make APTS working solution of desired concentration.

EXAMPLE DATA ANALYSIS AND FIGURES

Figure 1. Chemical structure for APTS [8-Aminopyrene-1,3,6-trisulfonic acid, trisodium salt] *CAS 196504-57-1*

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.