

# FluoroQuest™ TSA/PSA Antifade Mounting Medium

## \*Optimized for Tyramide and Styramide Imaging\*

Catalog Number: 44890

Unit Size: 5 mL

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Liquid
Molecular Weight	N/A
Soluble In	N/A

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

FluoroQuest™ TSA/PSA Antifade Mounting Medium is the only mounting medium that is optimized for tyramide (TSA) and styramide (PSA)-based fluorescence imaging applications. The antifade mounting medium is a specialized solution used in fluorescence microscopy to preserve the fluorescent signal emitted by fluorophores of TSA and PSA probes. When using fluorescent probes to label cellular components, it's crucial to preserve the fluorescence and prevent fading during the imaging process. Fluorescence fading occurs due to photobleaching, where the fluorophores lose their ability to emit light when exposed to intense repeated excitation light. The antifade mounting medium is designed to minimize photobleaching and maintain the fluorescence of the labeled structures over an extended period. This medium contains a combination of a few components that help protect the fluorophores and stabilize the fluorescent signal, including antioxidants, antifading agents, buffering agents and viscosity modifiers. AAT Bioquest also offers a few other mounting media for the general fluorescence imaging applications. It's essential to choose the appropriate antifade mounting medium based on the specific experimental needs and the fluorophores used. After mounting the sample with antifade medium, it is typically covered with a glass coverslip to protect it and then imaged using fluorescence microscopy to visualize and analyze the labeled structures.