

## XFD594 Phalloidin

Catalog Number: 23158

Unit Size: 300 Tests

### Product Details

---

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

### Chemical Properties

---

Appearance	Liquid purple
Molecular Weight	~1600
Soluble In	DMSO

### Spectral Properties

---

Excitation Wavelength	590 nm
Emission Wavelength	618 nm

### Applications

---

XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). XFD594 phalloidin conjugate is chemically equivalent to Alexa Fluor® 594 phalloidin. This red fluorescent phalloidin conjugate selectively binds to F-actins. Used at nanomolar concentrations, phalloidin derivatives are convenient probes for labeling, identifying and quantitating F-actins in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments. Fluorescent phalloidin derivatives have been used as an important tool in the study of actin networks at high resolution. AAT Bioquest offers a variety of fluorescent phalloidin derivatives with different colors for multicolor imaging applications.