

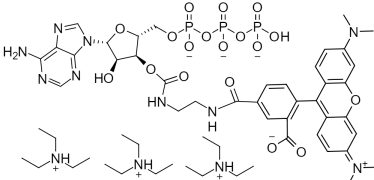
ATP-TAMRA conjugate [5-TAMRA-eda-ATP]

Catalog number: 13607
Unit size: 100 nmol

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	1128.13
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	552 nm
Emission Wavelength	578 nm

Applications

Fluorescently labeled ATP tracers are useful for monitoring ATP-binding proteins and other biomolecules. The ATP-TAMRA tracer was demonstrated to bind ATPases that are a class of enzymes that catalyze the decomposition of ATP into ADP and a free phosphate ion.