

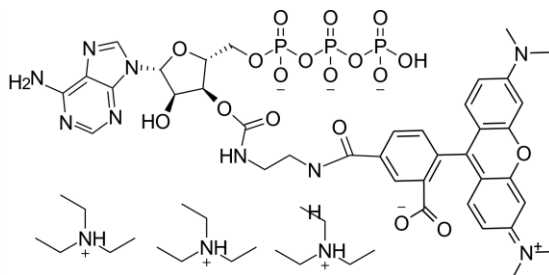
Product Information Sheet

Ordering Information

Product Number:	13607
Product Name:	ATP-TAMRA conjugate
Unit Size:	100 nmoles
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight:	1128.13
Appearance:	Red powder



Chemical Structure:	
Soluble in:	DMSO
Excitation Wavelength:	544 nm
Emission Wavelength:	575 nm

Application Notes

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.