

TSQ [N-(6-Methoxy-8-quinolyl)-p-toluenesulfonamide]Catalog number: 21254
Unit size: 5 mg**Product Details**

Storage Conditions Freeze (<math>-15\text{ }^{\circ}\text{C}</math>), Minimize light exposure

Expiration Date 12 months upon receiving

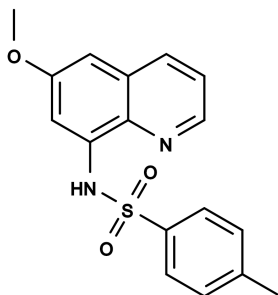
Chemical Properties

Appearance Off-white solid

Molecular Weight 328.38

Soluble In DMSO

Chemical Structure

**Applications**

TSQ is a fluorescent zinc indicator with high sensitivity.