

Tel: 408-733-1055; Fax: 408-733-1304 www.aatbio.com; E-Mail: info@aatbio.com

Product Information Sheet

Ordering Information

Product Number: 382

Product Name: 6-ROX [6-Carboxy-X-rhodamine] * Single isomers*

Size: 25 mg

Storage Conditions: <-15 °C and kept from light and moisture

Expiration Date: 12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 534.60

Appearance: Dark red solid

Chemical Structure:

Soluble in: DMSO
Excitation Wavelength: 570 nm
Emission Wavelength: 591 nm

Application Notes

6-ROX is the other purified single isomer of 5(6)-ROX. It is preferred for some complicated biological applications where reproducibility is more critical than material cost since the minor positional difference between 5-ROX and 6-ROX might significantly affect some biological properties of the underlying conjugates. 6-ROX is predominantly used to label nucleotides and for DNA sequencing. Currently AAT Bioquest is using a proprietary process to synthesize isomer-free 6-ROX without the conventional chromatographic purification. Cautions must be exercised to prevent the deterioration during storage.