

# Product Information

## Carboxy-X-Rhodamine, SE (ROX, SE)

Dye	Cat no.	Unit size	Ex/Em (nm)	MW
5(6)-ROX, SE	90005	25 mg	575/600	632
5-ROX, SE	90036	5 mg	568/595	632
6-ROX, SE	90037	5 mg	576/601	632
6-ROX, SE in DMSO (83 mg/mL, 0.13 M)	90098	100 uL	576/601	632
ROX-PEO8 SE, 5 isomer	90089	1 mg	568/595	1055

### Storage and Handling

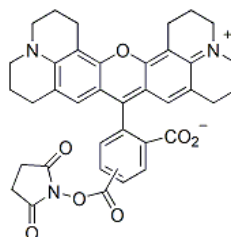
Store at -20°C desiccated and protect from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Stock solutions may be prepared in DMSO or DMF. Solutions can be aliquoted and stored with desiccant and protected from light at -20°C for at least a month.

### Product Description

Carboxy-X-rhodamine, SE (ROX, SE) is suitable for labeling oligonucleotides and is widely used in DNA sequencing and DNA probes based on Förster resonance energy transfer (FRET). For most applications, the 5- and 6-isomers of the dye can be used interchangeably. For some specialized applications, including Sanger sequencing, using a single isomer dye may be preferable to avoid altering the mobility of DNA fragments.

ROX-PEO8 SE, 5 isomer (Cat no. 90089) has a 32-angstrom water-soluble PEG spacer between the dye and the reactive group, to increase dye solubility and reduce potential steric interactions between the dye and the labeled molecule.

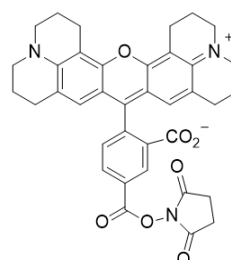


### 5(6)-ROX, SE

**Full chemical name:** 5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester

**Chemical Formula:** C<sub>37</sub>H<sub>33</sub>N<sub>3</sub>O<sub>7</sub>

**Color and Form:** Dark red solid

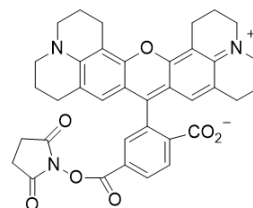


### 5-ROX, SE

**Full chemical name:** 5-Carboxy-X-rhodamine, succinimidyl ester, single isomer

**Chemical Formula:** C<sub>37</sub>H<sub>33</sub>N<sub>3</sub>O<sub>7</sub>

**Color and Form:** Dark red solid

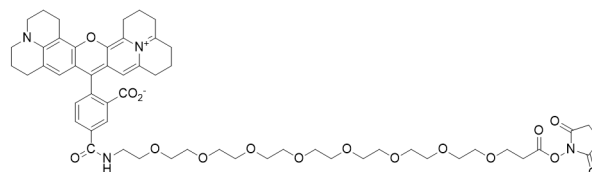


### 6-ROX, SE

**Full chemical name:** 6-Carboxy-X-rhodamine, succinimidyl ester, single isomer

**Chemical Formula:** C<sub>37</sub>H<sub>33</sub>N<sub>3</sub>O<sub>7</sub>

**Color and Form:** Dark red solid (Cat. no. 90037) or solution (Cat. no. 90098).



### 5-ROX-PEO8 SE

**Full chemical name:** 5-Carboxy-X-rhodamine-PEO<sub>8</sub>-propionate, succinimidyl ester

**Chemical Formula:** C<sub>56</sub>H<sub>70</sub>N<sub>4</sub>O<sub>16</sub>

**Color and Form:** Dark red solid

## Related Products

Catalog number	Product
90082	DMSO, Anhydrous
92103... 96078	CF® Dye Succinimidyl Ester
90016... 90049	Cyanine Dye Succinimidyl Ester
90017- 90019	Cyanine NS Dye Succinimidyl Ester
90034	5-TAMRA SE
90035	6-TAMRA SE
90097	6-TAMRA SE in DMSO
90022	5(6)-TAMRA SE (mixed isomers)
90081	Carboxyrhodamine 110-PEO4, SE, 5 isomer
90041	CFDA SE, 5 and 6-isomers
90020	FITC (Isomer I)
90029	FAM SE, 5 isomer
90023	FAM SE, 6 isomer
90024	FAM SE, 5 and 6 isomers
90085	FAM-PEO8 SE, 5 isomer
90018	TRITC, 5 isomer
90019	TRITC, 6 isomer
90017	TRITC, 5 and 6 isomers

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.